

Publication compilation

Compilation of articles performing BRET with Mithras or TriStar readers

Introduction

BRET (Bioluminescence Resonance Energy Transfer) is a powerful tool to study the interaction between two different molecules and is a popular method to study protein-protein interaction, especially in the field of G-protein coupled receptor research.

BRET requires the measurement of filtered luminescence, and benefits hence from microplate readers with high sensitivity in luminescence, such as the Mithras and TriStar series from Berthold Technologies. This document presents 180 articles published in peer-reviewed journals until June 2022, which use Mithras or TriStar readers to perform BRET, and include articles published in top journals such as Cell and Nature. The list is not exhaustive, but it provides many valuable references.

Year	Authors	Journal	Title	Link	Reader
2022	Chiara Borsari, Erhan Keles, Jacob A. McPhail, et al.	J. Am. Chem. Soc.	Covalent Proximity Scanning of a Distal Cysteine to Target PI3K α	https://doi.org/10.1021/jacs.1c13568	Mithras ² LB 943
2022	Farheen Badrealam Khan, Irfa Anwar, Elrashdy M. Redwan, et al.	J. of Dairy Science	Camel and bovine milk lactoferrins activate insulin receptor and its related AKT and ERK1/2 pathways	https://doi.org/10.3168/jds.2021-20934	TriStar ² LB 942
2022	Yeun Ju Kim, Nivedita Sengupta, Mira Sohn, et al.	EMBO Reports	Metabolic routing maintains the unique fatty acid composition of phosphoinositides	https://doi.org/10.15252/embr.202154532	TriStar ² LB 942
2022	Ana L. Moreno-Salinas, Brian J. Holleran, Estefania Y. Ojeda-Muñoz, et al.	Molecular Psychiatry	Convergent selective signaling impairment exposes the pathogenicity of latrophilin-3 missense variants linked to inheritable ADHD susceptibility	https://doi.org/10.1038/s41380-022-01537-3	TriStar ² LB 942
2022	Patrick Pagesy, Abdelouhab Bouaboud, Zhihao Feng, et al.	Cells	Short O-GlcNAcase Is Targeted to the Mitochondria and Regulates Mitochondrial Reactive Oxygen Species Level	https://doi.org/10.3390/cells11111827	TriStar ² LB 942
2022	Cécile Pétigny, Audrey-Ann Dumont, Hugo Giguère, et al.	Int. J. Mol. Sci.	Monitoring TRPC7 Conformational Changes by BRET Following GPCR Activation	https://doi.org/10.3390/ijms23052502	TriStar ² LB 942

2022	Jingjing Wang, Meng Wu, Zhangcheng Chen, et al.	Nat Commun.	The unconventional activation of the muscarinic acetylcholine receptor M4R by diverse ligands	https://doi.org/10.1038%2Fs41467-022-30595-y	Mithras LB 940
2021	Isra Al Zamel, Abdulrasheed Palakkott, Mohammed Akli Ayoub	FEBS Letters	Synergistic activation of thrombin and angiotensin II receptors revealed by bioluminescence resonance energy transfer	https://doi.org/10.1002/1873-3468.14187	TriStar ² LB 942
2021	Arshida Ashraf, Priti Mudgil, Abdulrasheed Palakkott, et al.	J. of Dairy Science	Molecular basis of the anti- diabetic properties of camel milk through profiling of its bioactive peptides on dipeptidyl peptidase IV (DPP-IV) and insulin receptor activity	https://doi.org/10.3168/jds.2020-18627	TriStar ² LB 942
2021	Bas Brouwers, Edson Mendes de Oliveira, Maria Martí-Solano, et al.	Cell Reports	Human MC4R variants affect endocytosis, trafficking and dimerization revealing multiple cellular mechanisms involved in weight regulation	https://doi.org/10.1016/j.celrep.2021.108862	Mithras LB 940
2021	Nidhi Chahal, Anjali Geethadevi, Surleen Kaura, et al.	Metabolism	Direct impact of gonadotropins on glucose uptake and storage in preovulatory granulosa cells: Implications in the pathogenesis of polycystic ovary syndrome	https://doi.org/10.1016/j.metabol.2020.154458	Mithras ² LB 943
2021	Xiaojing Cong, Damien Maurel, Hélène Déméné, et al.	Molecular Cell	Molecular insights into the biased signaling mechanism of the μ-opioid receptor	https://doi.org/10.1016/j.molcel.2021.07.033	Mithras LB 940
2021	J. Patrick Connick, James R. Reed, George F. Cawley, Wayne L. Backes	J. Biol. Chem.	Heme oxygenase-1 affects cytochrome P450 function through the formation of heteromeric complexes: Interactions between CYP1A2 and heme oxygenase-1	https://doi.org/10.1074/jbc.RA120.015911	TriStar LB 941
2021	Carlos Costas-Insua, Estefanía Moreno, Irene B. Maroto, et al.	J. Neurosci.	Identification of BiP as a CB1 Receptor-Interacting Protein That Fine-Tunes Cannabinoid Signaling in the Mouse Brain	https://dx.doi.org/10.1523%2FJNEUROSCI.0821-21.2021	Mithras LB 940
2021	Yann Chappe, Pauline Michel, Alexandre Joushomme, et al.	Molecular Pharmacology	High-Throughput Screening of TRPV1 Ligands in the Light of the Bioluminescence Resonance Energy Transfer Technique	https://doi.org/10.1124/molpharm.121.000271	TriStar ² LB 942
2021	Francesco De Pascali, Mohammed Akli Ayoub, Riccardo Benevelli, et al.	Int. J. Mol. Sci.	Pharmacological Characterization of Low Molecular Weight Biased Agonists at the Follicle Stimulating Hormone Receptor	https://dx.doi.org/10.3390%2Fijms22189850	Mithras ² LB 943
2021	Aliza T. Ehrlich, Meriem Semache, Pierre Couvineau, et al.	Mol. Brain	Ackr3-Venus knock-in mouse lights up brain vasculature	https://dx.doi.org/10.1186%2Fs13041-021-00862-y	Mithras LB 940, TriStar ² LB 942
2021	Chayma El Khamli, Laetitia Cobret, Jean-Michel Arrang, et al.	Int. J. Mol. Sci.	BRET Analysis of GPCR Dimers in Neurons and Non-Neuronal Cells: Evidence for Inactive, Agonist, and Constitutive Conformations	https://dx.doi.org/10.3390%2Fijms221910638	Mithras LB 940
2021	Nae Fu, Kazunori Sugiura, Kumiko Kondo, et al.	J. Biol. Chem.	Monitoring cellular redox dynamics using newly developed BRET-based redox sensor proteins	https://dx.doi.org/10.1016%2Fj.jbc.2021.101186	TriStar LB 941

2021	Alejandro Lillo, Jaume Lillo, Iu Raich, Cristina Miralpeix, et al.	Front. Cell. Neurosci.	Ghrelin and Cannabinoid Functional Interactions Mediated by Ghrelin/CB1 Receptor Heteromers That Are Upregulated in the Striatum From Offspring of Mice Under a High-Fat Diet	https://dx.doi.org/10.3389%2Ffncel.2021.786597	Mithras LB 940
2021	Jaume Lillo, Alejandro Lillo, David A. Zafra, et al.	Int. J. Mol. Sci.	Identification of the Ghrelin and Cannabinoid CB2 Receptor Heteromer Functionality and Marked Upregulation in Striatal Neurons from Offspring of Mice under a High-Fat Diet	https://dx.doi.org/10.3390%2Fijms22168928	Mithras LB 940
2021	Jing-Yu Lin, Zhao Yang, Chan Yang, et al.	J. Biol. Chem.	An ionic lock and a hydrophobic zipper mediate the coupling between an insect pheromone receptor BmOR3 and downstream effectors	https://dx.doi.org/10.1016%2Fj.jbc.2021.101160	Mithras LB 940
2021	Lei Liu, Zhiran Fan, Xavier Rovira, et al.	eLife	Allosteric ligands control the activation of a class C GPCR heterodimer by acting at the transmembrane interface	https://dx.doi.org/10.7554%2FeLife.70188	Mithras LB 940
2021	Shaoyong Lu, Xinheng He, Zhao Yang, et al.	Nat. Commun.	Activation pathway of a G protein-coupled receptor uncovers conformational intermediates as targets for allosteric drug design	https://dx.doi.org/10.1038%2Fs41467-021-25020-9	Mithras LB 940
2021	Xiaoyuan Ma, Eléonore W. E. Verweij, Marco Siderius, Rob Leurs, and Henry F. Vischer	Biomolecules	Identification of TSPAN4 as Novel Histamine H4 Receptor Interactor	https://dx.doi.org/10.3390%2Fbiom11081127	Mithras LB 940
2021	Mathilde Munier, Mohammed Ayoub, Valentine Suteau, et al.	Arch. Toxicol.	In vitro effects of the endocrine disruptor p,p'DDT on human choriogonadotropin/luteinizing hormone receptor signalling	https://doi.org/10.1007/s00204-021-03007-1	Mithras ² LB 943
2021	Wissem Deraredj Nadim, Shalina Hassanaly, Hélène Bénédicti, et al.	Biochem. Biophys. Res. Commun	The GTPase-activating protein-related domain of neurofibromin interacts with MC1R and regulates pigmentation-mediated signaling in human melanocytes	https://doi.org/10.1016/j.bbrc.2020.11.003	Mithras LB 940
2021	Shungo Nakamura, Nae Fu, Kumiko Kondo	J. Biol. Chem.	A luminescent Nanoluc-GFP fusion protein enables readout of cellular pH in photosynthetic organisms	https://doi.org/10.1074/jbc.RA120.016847	TriStar LB 941
2021	Yu-Qi Ping, Chunyou Mao, Peng Xiao, et al.	Nature	Structures of the glucocorticoid-bound adhesion receptor GPR97–Go complex	https://doi.org/10.1038/s41586-020-03083-w	Mithras LB 940
2021	Aratrika Saha, Patrick Connick, James R. Reed, et al.	Biochem. J.	Identification of the contact region responsible for the formation of the homomeric CYP1A2•CYP1A2 complex	https://doi.org/10.1042/BCJ20210269	TriStar LB 941
2021	Dexiu Wang, Yuliang Wang, Meiyang Shan, et al.	Experimental Cell Res.	Apelin receptor homodimer inhibits apoptosis in vascular dementia	https://doi.org/10.1016/j.yexcr.2021.112739	TriStar LB 941
2021	Shane C. Wright, Viktoriya Lukashova, Christian Le Gouill, et al.	PNAS	BRET-based effector membrane translocation assay monitors GPCR-promoted and endocytosis-mediated Gq activation at early endosomes	https://doi.org/10.1073/pnas.2025846118	TriStar ² LB 942

2021	Peng Xiao, Wei Yan, Lu Gou, et al.	Cell	Ligand recognition and allosteric regulation of DRD1-Gs signaling complexes	https://doi.org/10.1016/j.cell.2021.01.028	Mithras LB 940
2021	Paweł Zajdel, Katarzyna Grychowska, Szczepan Mogilski, et al.	J. Med. Chem.	Structure-Based Design and Optimization of FPPQ, a Dual-Acting 5-HT3 and 5-HT6 Receptor Antagonist with Antipsychotic and Procognitive Properties	https://dx.doi.org/10.1021%2Facs.jmedchem.1c00224	Mithras LB 940
2020	Charlotte Avet, Claudio Sturino, Sébastien Grastilleur, et al.	Communications Biology	The PAR2 inhibitor I-287 selectively targets Gαq and Gα12/13 signaling and has anti-inflammatory effects	https://doi.org/10.1038/s42003-020-01453-8	TriStar ² LB 942
2020	Xin Cai, Huannan Wang, Maochang Wang, et al.	Cellular Signaling	A novel phosphorylation site on orexin receptor 1 regulating orexinA-induced GRK2-biased signaling	https://doi.org/10.1016/j.cellsig.2020.109743	TriStar LB 941
2020	Valeria Francesconi, Elena Cichero, Evgeny V. Kanov, et al.	Pharmaceuticals	Novel 1-Amidino-4-Phenylpiperazines as Potent Agonists at Human TAAR1 Receptor: Rational Design, Synthesis, Biological Evaluation and Molecular Docking Studies	https://doi.org/10.3390/ph13110391	Mithras ² LB 943
2020	Thorsten Gnad, Gemma Navarro, Minna Lahesmaa, et al.	Cell Metab.	Adenosine/A2B Receptor Signaling Ameliorates the Effects of Aging and Counteracts Obesity	https://doi.org/10.1016/j.cmet.2020.06.006	Mithras LB 940
2020	Gergő Gulyás, Mira Sohn, Yeun Ju Kim, et al.	J Cell Sci	ORP3 phosphorylation regulates phosphatidylinositol 4-phosphate and Ca ²⁺ dynamics at plasma membrane–ER contact sites	https://doi.org/10.1242/jcs.237388	TriStar ² LB 942
2020	Bingyuan Ji, Liyan Shang, Chunmei Wang, et al.	Cellular Signaling	Roles for heterodimerization of APJ and B2R in promoting cell proliferation via ERK1/2-eNOS signaling pathway	https://doi.org/10.1016/j.cellsig.2020.109671	TriStar LB 941
2020	Hyunook Kang, Han-Sol Yang, Ah Young Ki, et al.	Structure	Conformational Dynamics and Functional Implications of Phosphorylated β-Arrestins	https://doi.org/10.1016/j.str.2019.12.008	TriStar ² LB 942
2020	Kuglae Kim, Tao Che, Ouliana Panova, et al.	Cell	Structure of a Hallucinogen-Activated Gq-Coupled 5-HT2A Serotonin Receptor	https://doi.org/10.1016/j.cell.2020.08.024	Mithras LB 940
2020	Thibaut Laboute, Jorge Gandía, Lucie P Pellissier, et al.	eLife	The orphan receptor GPR88 blunts the signaling of opioid receptors and multiple striatal GPCRs	https://doi.org/10.7554/eLife.50519	Mithras ² LB 943
2020	Viktoriya Lukasheva, Dominic Devost, Christian Le Gouill, et al.	Scientific Reports	Signal profiling of the β1AR reveals coupling to novel signalling pathways and distinct phenotypic responses mediated by β1AR and β2AR	https://doi.org/10.1038/s41598-020-65636-3	TriStar ² LB 942
2020	Maria Marti-Solano, Stephanie E. Crilly, Duccio Malinverni, et al.	Nature	Combinatorial expression of GPCR isoforms affects signalling and drug responses	https://doi.org/10.1038/s41586-020-2888-2	Mithras LB 940

2020	Tony Ngo, Bryan S. Stephens, Martin Gustavsson, et al.	PLOS Biology	Crosslinking-guided geometry of a complete CXCR2 receptor-chemokine complex and the basis of chemokine subfamily selectivity	https://doi.org/10.1371/journal.pbio.3000656	TriStar LB 941
2020	Sami Nourreddine, Geneviève Lavoie, Justine Paradis, et al.	Cell Reports	NF45 and NF90 Regulate Mitotic Gene Expression by Competing with Staufen-Mediated mRNA Decay	https://doi.org/10.1016/j.celrep.2020.107660	Mithras LB 940
2020	Joshua G. Pemberton, Yeun Ju Kim, Jana Humpolickova, et al.	J Cell Biol	Defining the subcellular distribution and metabolic channeling of phosphatidylinositol	https://doi.org/10.1083/jcb.201906130	TriStar ² LB 942
2020	Emmanuelle Poque, Delia Arnaud-Cormos, Lorenza Patrignoni, et al.	International Journal of Radiation Biology	Effects of radiofrequency fields on RAS and ERK kinases activity in live cells using the bioluminescence resonance energy transfer technique	https://doi.org/10.1080/09553002.2020.1730016	TriStar ² LB 942
2020	Lauren M. Slosky, Yushi Bai, Krisztian Toth, et al.	Cell	β -Arrestin-Biased Allosteric Modulator of NTSR1 Selectively Attenuates Addictive Behaviors	https://doi.org/10.1016/j.cell.2020.04.053	Mithras LB 940
2020	Reed M. Stein, Hye Jin Kang, John D. McCorvy,	Nature	Virtual discovery of melatonin receptor ligands to modulate circadian rhythms	https://doi.org/10.1038/s41586-020-2027-0	Mithras LB 940
2020	Fan Yang, Chunyou Mao, Lulu Guo, et al.	Nature	Structural basis of GPBAR activation and bile acid recognition	https://doi.org/10.1038/s41586-020-2569-1	Mithras LB 940
2019	Amanat Ali, Abdurashheed Palakkott, Arshida Ashraf, et al.	Front. Pharmacol.	Positive Modulation of Angiotensin II Type 1 Receptor-Mediated Signaling by LVV-Hemorphin-7	https://doi.org/10.3389/fphar.2019.01258	TriStar ² LB 942
2019	Takashi Baba, Daniel J Toth, Nivedita Sengupta, et al.	The EMBO Journal	Phosphatidylinositol 4,5-bisphosphate controls Rab7 and PLEKHM1 membrane cycling during autophagosome-lysosome fusion	https://doi.org/10.15252/embj.2018100312	TriStar ² LB 942
2019	Daniele Bolognini, Natasja Barki, Adrian J. Butcher, et al.	Nature Chemical Biology	Chemogenetics defines receptor-mediated functions of short chain free fatty acids	https://doi.org/10.1038/s41589-019-0270-1	TriStar LB 941
2019	Alexandre Connolly, Brian J. Holleran, Élie Simard, et al.	Biochemical Pharmacology	Interplay between intracellular loop 1 and helix VIII of the angiotensin II type 2 receptor controls its activation.	https://doi.org/10.1016/j.bcp.2019.07.018	TriStar ² LB 942
2019	Dany Fillion, Dominic Devost, Rory Sleno	Front. Endocrinol.	Asymmetric Recruitment of β -Arrestin1/2 by the Angiotensin II Type I and Prostaglandin F ₂ α Receptor Dimer	https://doi.org/10.3389/fendo.2019.00162	TriStar ² LB 942
2019	T. Chase Francis, Hideaki Yano, Tyler G. Demarest, et al.	Neuron	High-Frequency Activation of Nucleus Accumbens D1-MSNs Drives Excitatory Potentiation on D2-MSNs	https://doi.org/10.1016/j.neuron.2019.05.031	Mithras LB 940
2019	Jiankun Lyu, Sheng Wang, Trent E. Balius, et al.	Nature	Ultra-large library docking for discovering new chemotypes	https://doi.org/10.1038/s41586-019-0917-9	Mithras LB 940
2019	Nanako Nomura, Ryo Nishihara, Takahiro Nakajima, et al.	Anal. Chem.	Biothiol-Activatable Bioluminescent Coelenterazine Derivative for Molecular Imaging in Vitro and in Vivo	https://doi.org/10.1021/acs.analchem.9b00694	TriStar ² LB 942

2019	Louis-Philippe Picard, Anne-Marie Schonegge, and Michel Bouvier*	ACS Pharmacol. Transl. Sci.	Structural Insight into G Protein-Coupled Receptor Signaling Efficacy and Bias between Gs and β -Arrestin	https://doi.org/10.1021/acspstsci.9b00012	TriStar ² LB 942
2019	Mcheik Saria, Van Eeckhout Nils, De Poorter Cédric, et al.	Front. Immunol.	Coexpression of CCR7 and CXCR4 During B Cell Development Controls CXCR4 Responsiveness and Bone Marrow Homing	https://doi.org/10.3389/fimmu.2019.02970	Mithras LB 940
2019	Milan Sencanski, Sanja Glisic, Marko Šnajder, et al.	Scientific Reports	Computational design and characterization of nanobody-derived peptides that stabilize the active conformation of the β 2-adrenergic receptor (β 2-AR)	https://doi.org/10.1038/s41598-019-52934-8	TriStar ² LB 942
2019	Nivedita Sengupta Marko Jović Elena Barnaeva, et al.	Journal of Lipid Research	A large scale high-throughput screen identifies chemical inhibitors of phosphatidylinositol 4-kinase type II alpha[S]	https://doi.org/10.1194/jlr.D090159	TriStar ² LB 942
2019	Maja Susec, Milan Sencanski, Sanja Glisic, et al.	Neuropharmacology	Functional characterization of β 2-adrenergic and insulin receptor heteromers	https://doi.org/10.1016/j.neuropharm.2019.01.025	TriStar ² LB 942
2019	Qin-Qin Wang, Chun-Mei Wang, Bao-Hua Cheng, et al.	Cell Signaling	Signaling transduction regulated by 5-hydroxytryptamine 1A receptor and orexin receptor 2 heterodimers	https://doi.org/10.1016/j.cellsig.2018.11.014	TriStar LB 941
2018	Aleksander A. Aleksandrov, Veronika M. Knyazeva, Anna B. Volnova, et al.	Neurotoxicity Research	Identification of TAAR5 Agonist Activity of Alpha-NETA and Its Effect on Mismatch Negativity Amplitude in Awake Rats	https://doi.org/10.1007/s12640-018-9902-6	Mithras ² LB 943
2018	Vladimir Bobkov, Aurelien M. Zarca, Anneleen Van Hout, et al.	Biochemical Pharmacology	Nanobody-Fc constructs targeting chemokine receptor CXCR4 potently inhibit signaling and CXCR4-mediated HIV-entry and induce antibody effector functions	https://doi.org/10.1016/j.bcp.2018.10.014	Mithras LB 940
2018	J. Patrick Connick, James R. Reed and Wayne L. Backes	Drug Metabolism and Disposition	Characterization of Interactions Among CYP1A2, CYP2B4, and NADPH-cytochrome P450 Reductase: Identification of Specific Protein Complexes	https://doi.org/10.1124/dmd.117.078642	TriStar LB 941
2018	David Cotnoir-White, Mohamed El Ezzy, Pierre-Luc Boulay, et al.	PNAS	Monitoring ligand-dependent assembly of receptor ternary complexes in live cells by BRETfect	https://doi.org/10.1073/pnas.1716224115	TriStar ² LB 942
2018	Marie-Céline Frantz, Lucie P. Pellissier, Elsa Pflimlin, et al.	J. Med. Chem.	LIT-001, the First Nonpeptide Oxytocin Receptor Agonist that Improves Social Interaction in a Mouse Model of Autism	https://doi.org/10.1021/acs.jmedchem.8b00697	Mithras ² LB 943
2018	Martine Gonneau, Thierry Desprez, Marjolaine Martin, et al.	Current Biol.	Receptor Kinase THESEUS1 Is a Rapid Alkalinization Factor 34 Receptor in Arabidopsis	https://doi.org/10.1016/j.cub.2018.05.075	TriStar LB 941
2018	Manuel Grundmann, Nicole Merten, Davide Malfacini, et al.	Nature Commun.	Lack of beta-arrestin signaling in the absence of active G proteins	https://doi.org/10.1038/s41467-017-02661-3	Mithras ² LB 943
2018	Ryan D. Martin, Yalin Sun, Kyla Bourque, et al.	Cellular Signaling	Receptor- and cellular compartment-specific activation of the cAMP/PKA pathway by α 1-adrenergic and ETA endothelin receptors	https://doi.org/10.1016/j.cellsig.2018.01.002	TriStar

2018	Matthieu Masureel, Yaozhong Zou, Louis-Philippe Picard, et al.	Nature Chemical Biology	Structural insights into binding specificity, efficacy and bias of a β 2AR partial agonist	https://doi.org/10.1038/s41589-018-0145-x	TriStar ² LB 942
2018	Mira Sohn, Marek Korzeniowski, James P. Zewe, et al.	J Cell Biol	PI(4,5)P ₂ controls plasma membrane PI4P and PS levels via ORP5/8 recruitment to ER-PM contact sites	https://doi.org/10.1083/jcb.201710095	TriStar ² LB 942
2018	Qingjie Xue, Bo Bai, Bingyuan Ji, et al.	Front. Mol. Neurosci.	Ghrelin Through GHSR1a and OX1R Heterodimers Reveals a Gas-cAMP-cAMP Response Element Binding Protein Signaling Pathway in Vitro	https://doi.org/10.3389/fnmol.2018.00245	TriStar LB 941
2017	Nathalie Clement, Nicolas Renault, Jean-Luc Guillaume, et al.	Br. J. Pharmacol.	Importance of the second extracellular loop for melatonin MT1 receptor function and absence of melatonin binding in GPR50	https://doi.org/10.1111/bph.14029	Mithras LB 940
2017	Caroline Corbel, Sara Sartini, Elisabetta Levati, et al.	SLAS Discov.	Screening for Protein-Protein Interaction Inhibitors Using a Bioluminescence Resonance Energy Transfer (BRET)-Based Assay in Yeast	https://doi.org/10.1177/2472555216689530	Mithras LB 940
2017	Thorsten Fritzius, Rostislav Turecek, Riad Seddik, et al.	J. Neurosci.	KCTD Hetero-oligomers Confer Unique Kinetic Properties on Hippocampal GABAB Receptor-Induced K ⁺ Currents	https://doi.org/10.1523/JNEUROSCI.2181-16.2016	Mithras LB 940
2017	Arman Javadi, Ravi K Deevi, Emma Evergren, et al.	eLife	PTEN controls glandular morphogenesis through a juxtamembrane β -Arrestin1/ARHGAP21 scaffolding complex	https://doi.org/10.7554%2FeLife.24578	Mithras ² LB 943
2017	Bingyuan Ji, Haiqing Liu, Rumin Zhang, et al.	Cellular Signalling	Novel signaling of dynorphin at κ -opioid receptor/bradykinin B2 receptor heterodimers.	https://doi.org/10.1016/j.cellsig.2017.01.005	TriStar LB 941
2017	Laura Riccetti, Romain Yvinec, Danièle Klett, et al.	Sci. Rep.	Human Luteinizing Hormone and Chorionic Gonadotropin Display Biased Agonism at the LH and LH/CG Receptors	https://doi.org/10.1038/s41598-017-01078-8	Mithras ² LB 943
2017	Laura Riccetti, Danièle Klett, Mohammed Akli Ayoub, et al.	Molecular Human Reproduction	Heterogeneous hCG and hMG commercial preparations result in different intracellular signalling but induce a similar long-term progesterone response in vitro	https://doi.org/10.1093/molehr/gax047	Mithras ² LB 943
2017	William G. Robichaux III, Melissa Branham-O'Connor, Il-Young Hwang, et al.	Journal of Pharmacology and Experimental Therapeutics	Regulation of Chemokine Signal Integration by Activator of G-Protein Signaling 4 (AGS4)	https://doi.org/10.1124/jpet.116.238436	TriStar LB 941
2017	Hermanus Johannes Ruigrok, Guillaume Shahid, Bertrand Goudeau, et al.	Biophys. J.	Full-Spectral Multiplexing of Bioluminescence Resonance Energy Transfer in Three TRPV Channels	https://doi.org/10.1016/j.bpj.2016.11.3197	TriStar ² LB 942
2017	Thomas H. Smith, Julia G. Li, Michael R. Dore, JoAnn Trejo	J. Biol. Chem.	Protease-activated receptor-4 and purinergic receptor P2Y ₁₂ dimerize, co-internalize, and activate Akt signaling via endosomal recruitment of β -arrestin	https://doi.org/10.1074/jbc.M117.782359	TriStar LB 941

2016	Abdulrasheed O. Abdulrahman, Mohammad A. Ismael, Khaled Al-Hosaini, et al.	Front. Endocrinol.	Differential Effects of Camel Milk on Insulin Receptor Signaling – Toward Understanding the Insulin-Like Properties of Camel Milk	https://doi.org/10.3389/fendo.2016.00004	Mithras ² LB 943
2016	Mohammed Akli Ayoub, Romain Yvinec, Gwenhaël Jégota, et al.	Mol. Cell. Endocrinol.	Profiling of FSHR negative allosteric modulators on LH/CGR reveals biased antagonism with implications in steroidogenesis	https://doi.org/10.1016/j.mce.2016.07.013	Mithras ² LB 943
2016	Constance Auvynet, Camille Baudesson de Chanville, Patricia Hermand, et al.	FASEB J.	ECL1i, d(LGTLKLC), a novel, small peptide that specifically inhibits CCL2-dependent migration	https://doi.org/10.1096/fj.201500116	TriStar
2016	Cédric M. Blouin, Yannick Hamon, Pauline Gonnord, et al.	Cell	Glycosylation-Dependent IFN- γ R Partitioning in Lipid and Actin Nanodomains Is Critical for JAK Activation	https://doi.org/10.1016/j.cell.2016.07.003	Mithras LB 940
2016	Sophie J. Bradley, Coen H. Wiegman, Max Maza Iglesias, et al.	PNAS	Mapping physiological G protein-coupled receptor signaling pathways reveals a role for receptor phosphorylation in airway contraction	https://doi.org/10.1073/pnas.1521706113	TriStar LB 941
2016	Nicole E. Brown, Nevin A. Lambert, John R. Hepler	Pharmacology Research & Perspectives	RGS14 regulates the lifetime of G α -GTP signaling but does not prolong G $\beta\gamma$ signaling following receptor activation in live cells	https://doi.org/10.1002/prp2.249	TriStar LB 941
2016	Nicolas Boute, Peter Lowe, Sven Berger, et al.	Front. Pharmacol.	NanoLuc Luciferase – A Multifunctional Tool for High Throughput Antibody Screening	https://doi.org/10.3389/fphar.2016.00027	Mithras LB 940
2016	Xin Cai, Bo Bai, Rumin Zhang, et al.	Sci. Rep.	Apelin receptor homodimer-oligomers revealed by single-molecule imaging and novel G protein-dependent signaling	https://doi.org/10.1038/srep40335	TriStar LB 941
2016	Alexandra Gonzalez, Matthieu Broussas, Charlotte Beau-Larvor, et al.	Int. J. Cancer	A novel antagonist anti-cMet antibody with antitumor activities targeting both ligand-dependent and ligand-independent c-Met receptors	https://doi.org/10.1002/ijc.30174	Mithras LB 940
2016	Elise Goyet, Nathalie Bouquier, Vincent Ollendorff & Julie Perroy	Sci. Rep.	Fast and high resolution single-cell BRET imaging.	https://doi.org/10.1038/srep28231	Mithras LB 940
2016	Valentina Kubale, Kaja Blagotinšek, Jane Nøhr, et al.	Int. J. Mol. Sci.	The Conserved Arginine Cluster in the Insert of the Third Cytoplasmic Loop of the Long Form of the D2 Dopamine Receptor (D2L-R) Acts as an Intracellular Retention Signal	https://doi.org/10.3390/ijms17071152	TriStar ² LB 942
2016	Mi-Hye Lee, Kathryn M. Appleton, Erik G. Strungs, et al.	Nature	The conformational signature of β -arrestin2 predicts its trafficking and signalling functions	https://doi.org/10.1038/nature17154	TriStar LB 941
2016	Aashish Manglik, Henry Lin, Dipendra K. Aryal, et al.	Nature	Structure-based discovery of opioid analgesics with reduced side effects	https://doi.org/10.1038/nature19112	Mithras LB 940
2016	Wisse Deraredj Nadim, Séverine Chaumont-Dubel, Fahima Madouri, et al.	PNAS	Physical interaction between neurofibromin and serotonin 5-HT ₆ receptor promotes receptor constitutive activity	https://doi.org/10.1073/pnas.1600914113	Mithras LB 940

2016	Sarah M. Pedersen, Waipan Chan, Rakhi P. Jattani, et al.	Molecular and Cellular Biology	Negative Regulation of CARD11 Signaling and Lymphoma Cell Survival by the E3 Ubiquitin Ligase RNF181	https://doi.org/10.1128/MCB.00876-15	TriStar LB 941
2016	Amynah A. Pradhan, Julie Perroy, Wendy M. Walwyn, et al.	J. Neurosci.	Agonist-Specific Recruitment of Arrestin Isoforms Differentially Modify Delta Opioid Receptor Function	https://doi.org/10.1523/JNEUROSCI.4124-15.2016	Mithras LB 940
2016	Mira Sohn, Pavlina Ivanova, H. Alex Brown, et al.	PNAS	Lenz-Majewski mutations in PTDSS1 affect phosphatidylinositol 4-phosphate metabolism at ER-PM and ER-Golgi junctions	https://doi.org/10.1073/pnas.1525719113	TriStar ² LB 942
2015	Mohammed Akli Ayoub, Flavie Landomiel, Nathalie Gallay, et al.	Front. Endocrinol.	Assessing gonadotropin receptor function by resonance energy transfer-based assays	https://doi.org/10.3389/fendo.2015.00130	Mithras ² LB 943
2015	Nicole E. Brown, Devrishi Goswami, Mary Rose Branch, et al.	J. Biol. Chem.	Integration of G Protein α (G α) Signaling by the Regulator of G Protein Signaling 14 (RGS14)	https://doi.org/10.1074/jbc.M114.634329	TriStar LB 941
2015	Buxin Chen, Antonio G. Soto, Luisa J. Coronel, et al.	Molecular Pharmacology	Characterization of Thrombin-Bound Dabigatran Effects on Protease-Activated Receptor-1 Expression and Signaling In Vitro	https://doi.org/10.1124/mol.114.096446	TriStar LB 941
2015	Julie Davies, Jing Chen, Ryan Pink, et al.	Sci. Rep.	Orexin receptors exert a neuroprotective effect in Alzheimer's disease (AD) via heterodimerization with GPR103.	https://doi.org/10.1038/srep12584	TriStar LB 941
2015	Michael R. Does, Huilan Lin, Neil J. Grimsey, et al.	Molecular Biology of the Cell	The α -arrestin ARDC3 mediates ALIX ubiquitination and G protein-coupled receptor lysosomal sorting	https://doi.org/10.1091/mbc.E15-05-0284	TriStar LB 941
2015	Tien-Hung Lan, Qiuju Liu, Chunman Li, et al.	Sci. Rep.	BRET evidence that β 2 adrenergic receptors do not oligomerize in cells	https://doi.org/10.1038/srep10166	Mithras LB 940
2015	Rodrigo Mendoza-Sanchez, David Cotnoir-White, Justyna Kulpa, et al.	Bioorganic & Medicinal Chemistry	Design, synthesis and evaluation of antiestrogen and histone deacetylase inhibitor molecular hybrids	https://doi.org/10.1016/j.bmc.2015.11.005	TriStar LB 941
2015	Yuki Ohba, Michio Nakaya, Kenji Watari, et al.	Biochem. Biophys. Res. Commun.	GRK6 phosphorylates I κ B α at Ser ³² /Ser ³⁶ and enhances TNF- α -induced inflammation	https://doi.org/10.1016/j.bbrc.2015.04.027	Mithras LB 940
2015	William G. Robichaux III, Sukru S. Oner, Stephen M. Lanier and Joe B. Blumer	Molecular Pharmacology	Direct Coupling of a Seven-Transmembrane-Span Receptor to a G α i G-Protein Regulatory Motif Complex	https://doi.org/10.1124/mol.115.097741	TriStar LB 941
2015	Ramona Schrage, Anna-Lena Schmitz, Evelyn Gaffal, et al.	Nature Commun.	The experimental power of FR900359 to study Gq-regulated biological processes	https://doi.org/10.1038/ncomms10156	Mithras LB 940
2014	Sylvain Armando, Julie Quoyer, Viktorya Lukashova, et al.	The FASEB Journal	The chemokine CXCL4 and CXCL2 receptors form homo- and heterooligomers that can engage their signaling G-protein effectors and β arrestin.	https://doi.org/10.1096/fj.13-242446	Mithras LB 940, TriStar ² LB 942

2014	Bo Bai, Xin Cai, Yunlu Jiang, et al.	Journal of Cellular and Molecular Medicine	Heterodimerization of apelin receptor and neurotensin receptor 1 induces phosphorylation of ERK1/2 and cell proliferation via Gαq-mediated mechanism	https://doi.org/10.1111/jcmm.12404	TriStar LB 941
2014	Marianne Bjordal, Nathalie Arquier, Julie Kniazeff, et al.	Cell	Sensing of Amino Acids in a Dopaminergic Circuitry Promotes Rejection of an Incomplete Diet in Drosophila	https://doi.org/10.1016/j.cell.2013.12.024	Mithras LB 940
2014	Ghil, Sungho, Sungho Ghil, Kelly L. McCoy, John R. Hepler	PloS one	Regulator of G protein signaling 2 (RGS2) and RGS4 form distinct G protein-dependent complexes with protease activated-receptor 1 (PAR1) in live cells.	https://doi.org/10.1371/journal.pone.0095355	TriStar LB 941
2014	Naïma Hanoun, Samuel Fritsch, Odile Gayet, et al.	J. Biol. Chem.	The E3 Ubiquitin Ligase Thyroid Hormone Receptor-interacting Protein 12 Targets Pancreas Transcription Factor 1a for Proteasomal Degradation	https://doi.org/10.1074/jbc.M114.620104	TriStar LB 941
2014	Gerda M. Hübner, Jane Nøhr Larsen, Barbara Guerra, et al.	Mol. Cell. Biochem.	Evidence for aggregation of protein kinase CK2 in the cell: a novel strategy for studying CK2 holoenzyme interaction by BRET2	https://doi.org/10.1007/s11010-014-2196-y	TriStar ² LB 942
2014	J-H Kang, S. A. Hassan, P. Zhao, et al.	Biochim. Biophys. Acta	Impact of subdomain D1 of the short form S1b of the human prolactin receptor on its inhibitory action on the function of the long form of the receptor induced by prolactin	https://dx.doi.org/10.1016%2Fj.bbagen.2014.04.006	Mithras LB 940
2014	Stefanie L. Makowski, Zhaoquan Wang, Joel L. Pomerantz	Molecular and Cellular Biology	A Protease-Independent Function for SPPL3 in NFAT Activation	https://doi.org/10.1128/MCB.01124-14	TriStar LB 941
2014	Anna-Lena Schmitz, Ramona Schrage, Evelyn Gaffal, et al.	Cell Chem. Biol.	A Cell-Permeable Inhibitor to Trap Gαq Proteins in the Empty Pocket Conformation	https://doi.org/10.1016/j.chembiol.2014.06.003	Mithras LB 940
2014	Salvador Sierra, Natasha Luquin, Alberto J. Rico, et al.	Brain Struct. Funct.	Detection of cannabinoid receptors CB1 and CB2 within basal ganglia output neurons in macaques: changes following experimental parkinsonism	https://doi.org/10.1007/s00429-014-0823-8	Mithras LB 940
2014	Yolanda Williams-Bey, Cedric Boullaran, Ali Vural, et al.	PLoS One	Omega-3 Free Fatty Acids Suppress Macrophage Inflammasome Activation by Inhibiting NF-κB Activation and Enhancing Autophagy	https://doi.org/10.1371/journal.pone.0097957	Mithras LB 940
2013	Huilan Lin, JoAnn Trejo	J. Biol. Chem.	Transactivation of the PAR1-PAR2 Heterodimer by Thrombin Elicits β-Arrestin-mediated Endosomal Signaling	https://doi.org/10.1074/jbc.M112.439950	TriStar LB 941
2013	Van B. Lu, Henry L. Puhl III and Stephen R. Ikeda	Molecular Pharmacology	N-Arachidonyl Glycine Does Not Activate G Protein–Coupled Receptor 18 Signaling via Canonical Pathways	https://doi.org/10.1124/mol.112.081182	TriStar LB 941
2013	Kasturi Pal, Maneesh Mathur, Puneet Kumar, Kathryn DeFea	J. Biol. Chem.	Divergent β-Arrestin-dependent Signaling Events Are Dependent upon Sequences within G-protein-coupled Receptor C Termini	https://doi.org/10.1074/jbc.M112.400234	TriStar LB 941

2013	Christopher P. Vellano, Nicole E. Brown, Joe B. Blumer, John R. Hepler	J. Biol. Chem.	Assembly and Function of the Regulator of G protein Signaling 14 (RGS14)-H-Ras Signaling Complex in Live Cells Are Regulated by Gαi1 and Gαi-linked G Protein-coupled Receptors	https://doi.org/10.1074/jbc.M112.440057	TriStar LB 941
2013	Yu-Jin Won, Van B. Lu, Henry L. Puhl III and Stephen R. Ikeda	J. Neurosci.	β-Hydroxybutyrate Modulates N-Type Calcium Channels in Rat Sympathetic Neurons by Acting as an Agonist for the G-Protein-Coupled Receptor FFA3	https://doi.org/10.1523/JNEUROSCI.3102-13.2013	TriStar LB 941
2012	Larry S. Barak and Sean Peterson	Biochemistry	Modeling of Bias for the Analysis of Receptor Signaling in Biochemical Systems	https://doi.org/10.1021/bi201308s	Mithras LB 940
2012	Andreas Bock, Nicole Merten, Ramona Schrage, et al.	Nat. Commun.	The allosteric vestibule of a seven transmembrane helical receptor controls G-protein coupling	https://doi.org/10.1038/ncomms2028	Mithras LB 940
2012	Nikolaj Kulahin, Samra J. Sanni, Rita Slaaby, et al.	Research J. Recept. Signal Transduct.	A BRET assay for monitoring insulin receptor interactions and ligand pharmacology	https://doi.org/10.3109/10799893.2011.647351	Mithras LB 940
2012	Tien-Hung Lan, Qiuju Liu, Chunman Li, et al.	Traffic	Sensitive and High Resolution Localization and Tracking of Membrane Proteins in Live Cells with BRET	https://doi.org/10.1111/j.1600-0854.2012.01401.x	Mithras LB 940
2012	R. Quintana-Portillo, A. Canfrán-Duque, T. Issad, et al.	Biochemical Pharmacology	Sam68 interacts with IRS1	https://doi.org/10.1016/j.bcp.2011.09.030	TriStar LB 941
2012	James R. Reed, J. Patrick Connick, Dongmei Cheng, et al.	Biochem. J.	Effect of homomeric P450–P450 complexes on P450 function	https://doi.org/10.1042/BJ20120636	TriStar LB 941
2012	M Seillier, S Peugeot, O Gayet, et al.	Cell Death & Differentiation	TP53INP1, a tumor suppressor, interacts with LC3 and ATG8-family proteins through the LC3-interacting region (LIR) and promotes autophagy-dependent cell death	https://doi.org/10.1038/cdd.2012.30	TriStar LB 941
2012	Hideo Takakura, Mitsuru Hattori, Masaki Takeuchi, and Takeaki Ozawa	ACS Chem. Biol.	Visualization and Quantitative Analysis of G Protein-Coupled Receptor–β-Arrestin Interaction in Single Cells and Specific Organs of Living Mice Using Split Luciferase Complementation	https://doi.org/10.1021/cb200360z	TriStar LB 941
2012	S. Vasseur, S. Afzal, R. Tomasini, et al.	Oncogene	Consequences of DJ-1 upregulation following p53 loss and cell transformation	https://doi.org/10.1038/onc.2011.268	TriStar LB 941
2011	Etha S. Burstein, Maria L. Carlsson, Michelle Owens, et al.	J. Neural Transm.	In vitro evidence that (–)-OSU6162 and (+)-OSU6162 produce their behavioral effects through 5-HT _{2A} serotonin and D ₂ dopamine receptors	https://doi.org/10.1007/s00702-011-0701-y	Mithras LB 940
2011	Caroline Corbel, Qian Wang, Hadjira Bousserouel, et al.	Biotechnol. J.	First BRET-based screening assay performed in budding yeast leads to the discovery of CDK5/p25 interaction inhibitors	https://doi.org/10.1002/biot.201100138	Mithras LB 940
2011	Christopher C. Marohnic, Warren J. Huber III, J. Patrick Connick	Archives of Biochemistry and Biophysics	Mutations of human cytochrome P450 reductase differentially modulate heme oxygenase-1 activity and oligomerization	https://doi.org/10.1016/j.abb.2011.06.008	TriStar LB 941

2011	Christopher P. Vellano, Ellen M. Maher, John R. Hepler, and Joe B. Blumer	J. Biol. Chem.	G Protein-coupled Receptors and Resistance to Inhibitors of Cholinesterase-8A (Ric-8A) Both Regulate the Regulator of G Protein Signaling 14 (RGS14)-Gαi1 Complex in Live Cells	https://doi.org/10.1074/jbc.M111.274928	TriStar LB 941
2011	Fallou Wade, Agathe Espagne, Marie-Annick Persuy, et al.	J. Biol. Chem.	Relationship between Homologous oligomerization of a Mammalian Olfactory Receptor and Its Activation State Demonstrated by Bioluminescence Resonance Energy Transfer	https://doi.org/10.1074/jbc.M110.184580	TriStar LB 941
2010	T Aoki, S Shimizu, E Urano, Y Futahashi, et al.	Gene Therapy	Improvement of lentiviral vector-mediated gene transduction by genetic engineering of the structural protein Pr55Gag	https://doi.org/10.1038/gt.2010.61	TriStar LB 941
2010	Johan Bacart, Audrey Leloire, Angélique Levoe, et al.	FEBS Lett.	Evidence for leptin receptor isoforms heteromerization at the cell surface	https://doi.org/10.1016/j.febslet.2010.03.033	Mithras LB 940
2010	Billy Breton, Étienne Sauvageau, Joris Zhou, et al.	Biophys. J.	Multiplexing of Multicolor Bioluminescence Resonance Energy Transfer	https://doi.org/10.1016/j.bpj.2010.10.025	Mithras LB 940
2010	Karima Bourougaa, Nadia Naski, Cedric Boularan, et al.	Molecular Cell	Endoplasmic Reticulum Stress Induces G2 Cell-Cycle Arrest via mRNA Translation of the p53 Isoform p53/47	https://doi.org/10.1016/j.molcel.2010.01.041	Mithras LB 940
2010	Shan Cen, Zong-Gen Peng, Xiao-Yu Li, et al.	J. Bio. Chem.	Small Molecular Compounds Inhibit HIV-1 Replication through Specifically Stabilizing APOBEC3G	https://doi.org/10.1074/jbc.M109.085308	Mithras LB 940
2010	Luc De Vries, Frédéric Finana, Frédéric Cachoux, et al.	Cellular Signaling	Cellular BRET assay suggests a conformational rearrangement of preformed TrkB/Shc complexes following BDNF-dependent activation	https://doi.org/10.1016/j.cellsig.2009.09.027	Mithras LB 940
2010	Stéphanie Gravel, Camille Malouf, Philip E. Boulais, et al.	J. Biol. Chem.	The Peptidomimetic CXCR4 Antagonist TC14012 Recruits β-Arrestin to CXCR7: ROLES OF RECEPTOR DOMAINS*	https://doi.org/10.1074/jbc.C110.147470	Mithras LB 940
2010	Sudhakaranmayi Kuravi, Tien-Hung Lan, Arnab Barik, et al.	Biophys. J.	Third-Party Bioluminescence Resonance Energy Transfer Indicates Constitutive Association of Membrane Proteins: Application to Class A G-Protein-Coupled Receptors and G-Proteins	https://doi.org/10.1016/j.bpj.2010.02.004	Mithras LB 940
2010	Anne Mazars, Robin Fähræus	Biotechnol. J.	Using BRET to study chemical compound-induced disruptions of the p53-HDM2 interactions in live cells	https://doi.org/10.1002/biot.200900272	Mithras LB 940
2010	Naomi Misawa, A. K. M. Kafi, Mitsuru Hattori, et al.	Anal. Chem.	Rapid and High-Sensitivity Cell-Based Assays of Protein-Protein Interactions Using Split Click Beetle Luciferase Complementation: An Approach to the Study of G-Protein-Coupled Receptors	https://doi.org/10.1021/ac100104g	TriStar LB 941
2010	Sukru Sadik Oner, Ningfei An, Ali Vural, et al.	J. Biol. Chem.	Regulation of the AGS3-Gαi Signaling Complex by a Seven-transmembrane Span Receptor	https://doi.org/10.1074/jbc.M110.138073	TriStar LB 941

2010	Sukru Sadik Oner, Ellen M. Maher, Billy Breton, et al.	J. Biol. Chem.	Receptor-regulated interaction of activator of G-protein signaling-4 and Gai	https://doi.org/10.1074/jbc.C109.088070	TriStar LB 941
2009	Nuria Cabello, Jorge Gandía, Daniela C. G. Bertarelli, et al.	J. Neurochem.	Metabotropic glutamate type 5, dopamine D2 and adenosine A2a receptors form higher-order oligomers in living cells	https://doi.org/10.1111/j.1471-4159.2009.06078.x	Mithras LB 940
2009	Makiko Hamatake, Toru Aoki, Yuko Futahashi, et al.	Cancer Sci.	Ligand-independent higher-order multimerization of CXCR4, a G-protein-coupled chemokine receptor involved in targeted metastasis	https://doi.org/10.1111/j.1349-7006.2008.00997.x	TriStar LB 941
2009	Maud Kamal, Marcel Marquez, Virginie Vauthier, et al.	Biotechnol. J.	Improved donor/acceptor BRET couples for monitoring β -arrestin recruitment to G protein-coupled receptors	https://doi.org/10.1002/biot.200900016	Mithras LB 940
2009	Angélique Levoe, Karl Balabanian, Françoise Baleux, et al.	Blood	CXCR7 heterodimerizes with CXCR4 and regulates CXCL12-mediated G protein signaling	https://doi.org/10.1182/blood-2008-12-196618	Mithras LB 940
2009	Marie-Laure Rives, Claire Vol, Yugo Fukazawa, et al.	EMBO J.	Crosstalk between GABAB and mGlu1a receptors reveals new insight into GPCR signal integration	https://doi.org/10.1038/emboj.2009.177	Mithras LB 940
2008	Larry S. Barak, Ali Salahpour, Xiaodong Zhang, et al.	Mol Pharmacol	Pharmacological Characterization of Membrane-Expressed Human Trace Amine-Associated Receptor 1 (TAAR1) by a Bioluminescence Resonance Energy Transfer cAMP Biosensor	https://dx.doi.org/10.1124%2Fmol.108.048884	Mithras LB 940
2008	Paulina Carriba, Gemma Navarro, Francisco Ciruela,	Nature Methods	Detection of heteromerization of more than two proteins by sequential BRET-FRET	https://doi.org/10.1038/nmeth.1229	Mithras LB 940
2008	Jorge Gandia, Jorge Galino, Olavo B. Amaral, et al.	FEBS Letters	Detection of higher-order G protein-coupled receptor oligomers by a combined BRET-BiFC technique	https://doi.org/10.1016/j.febslet.2008.07.045	Mithras LB 940
2008	Bernard Masri, Ali Salahpour, Michael Didriksen, et al.	PNAS	Antagonism of dopamine D2 receptor/ β -arrestin 2 interaction is a common property of clinically effective antipsychotics	https://doi.org/10.1073/pnas.0803522105	Mithras LB 940
2008	Gemma Navarro, Paulina Carriba, Jorge Gandí, et al.	Sci. World	Detection of Heteromers Formed by Cannabinoid CB1, Dopamine D2, and Adenosine A2A G-Protein-Coupled Receptors by Combining Bimolecular Fluorescence Complementation and Bioluminescence Energy Transfer	https://doi.org/10.1100/tsw.2008.136	Mithras LB 940
2008	Julie Perroy, Fabrice Raynaud, Vincent Homburger, et al.	J. Biol. Chem.	Direct Interaction Enables Cross-talk between Ionotropic and Group I Metabotropic Glutamate Receptors	https://doi.org/10.1074/jbc.M705661200	Mithras LB 940
2007	Mohammed A. Ayoub, Damien Maurel, Virginie Binet, et al.	Molecular Pharmacology	Real-Time Analysis of Agonist-Induced Activation of Protease-Activated Receptor 1/Gai1 Protein Complex Measured by Bioluminescence Resonance Energy Transfer in Living Cells	https://doi.org/10.1124/mol.106.030304	Mithras LB 940

2007	Randy L. Bogan, Tracy L. Davis, Gordon D. Niswender	J. Steroid Biochem. Mol. Biol	Peripheral-type benzodiazepine receptor (PBR) aggregation and absence of steroidogenic acute regulatory protein (StAR)/PBR association in the mitochondrial membrane as determined by bioluminescence resonance energy transfer (BRET)	https://doi.org/10.1016/j.jsbmb.2006.10.007	Mithras LB 940
2007	Cédric Boularan, Mark G. H. Scott, Karima Bourougaa, et al.	PNAS	β -arrestin 2 oligomerization controls the Mdm2-dependent inhibition of p53	https://doi.org/10.1073/pnas.0705550104	Mithras LB 940
2007	Cyril Couturier, Chamsy Sarkis, Karin Séron, et al.	PNAS	Silencing of OB-RGRP in mouse hypothalamic arcuate nucleus increases leptin receptor signaling and prevents diet- induced obesity	https://doi.org/10.1073/pnas.0706671104	Mithras LB 940
2007	Hans H Schiffer, Esther C Reding, Stephen R Fuhs, et al.	Mol. Pharmacol.	Pharmacology and signaling properties of epidermal growth factor receptor isoforms studied by bioluminescence resonance energy transfer	https://doi.org/10.1124/mol.106.027656	Mithras LB 940
2006	Fadi F. Hamdan, Yann Percherancier, Billy Breton and Michel Bouvier	Current Protocols in Neuroscience	Monitoring Protein-Protein Interactions in Living Cells by Bioluminescence Resonance Energy Transfer (BRET)	https://doi.org/10.1002/0471142301.ns0523s34	Mithras LB 940
2006	Kevin D. G. Pflieger, Jasmin R. Dromey, Matthew B. Dalrymple, et al.	Cellular Signaling	Extended bioluminescence resonance energy transfer (eBRET) for monitoring prolonged protein–protein interactions in live cells	https://doi.org/10.1016/j.cellsig.2006.01.004	Mithras LB 940
2005	Pascale G Charest, Sonia Terrillon and Michel Bouvier	EMBO Reports	Monitoring agonist-promoted conformational changes of β - arrestin in living cells by intramolecular BRET	https://doi.org/10.1038/sj.embor.7400373	Mithras LB 940
2005	Céline Galés, R Victor Rebois, Mireille Hogue, et al.	Nature Methods	Real-time monitoring of receptor and G-protein interactions in living cells	https://doi.org/10.1038/nmeth743	Mithras LB 940
2005	Evi Kostenis, Graeme Milligan, Arthur Christopoulos, et al.	Circulation	G-Protein–Coupled Receptor Mas Is a Physiological Antagonist of the Angiotensin II Type 1 Receptor	https://doi.org/10.1161/01.CIR.0000160867.23556.7D	Mithras LB 940
2005	Shirley Wilson, Graeme Wilkinson and Graeme Milligan	J. Biol. Chem.	The CXCR1 and CXCR2 Receptors Form Constitutive Homo- and Heterodimers Selectively and with Equal Apparent Affinities*	https://doi.org/10.1074/jbc.M413475200	Mithras LB 940
2004	Julie Perroy, Stephanie Pontier, Pascale G Charest, Muriel Aubry & Michel Bouvier	Nature Methods	Real-time monitoring of ubiquitination in living cells by BRET	https://doi.org/10.1038/nmeth722	Mithras LB 940

For Research Use Only. Not for use in diagnostic procedures.

© 2022 Berthold Technologies. All rights reserved. The trademarks mentioned herein are the property of Berthold Technologies or their respective owners.

Berthold Technologies GmbH & Co. KG. www.berthold.com