

Newsletter April 2010

## SBS & ISBC Issue

### Newsletter Topics

- ▶ **Brain Absorption Assay in ADME/Tox studies**
- ▶ **Berthold Technologies at SBS, ISBC & EUCARPIA**
- ▶ **Featured Product: NightSHADE Plant Imager**
- ▶ **Berthold on Facebook**
- ▶ **New Application & Technical Notes**
- ▶ **Recent Publications**

### Mithras plate reader for ADME/Tox assays

The [Transil® Brain Absorption Assay](#) from Sovicell has been successfully performed with the [Mithras](#) multimode reader. The kit is designed to assess drug candidates' ability to cross the blood brain barrier. The assay determines drug candidates' affinities to brain membranes and thus the brain free fraction available to interact with receptors.

These parameters may be used to predict whether a candidate has CNS drug like properties or whether it is likely to show CNS side effects.

### Berthold Technologies present at SBS and ISBC and EUCARPIA

At [SBS](#) 2010 conference of the Society for Biomolecular Sciences taking place in Phoenix, Arizona, Berthold Technologies will present its [Mithras](#) multimode and [CentroPRO](#) luminescence microplate readers.

Please come and see us at booth 603.

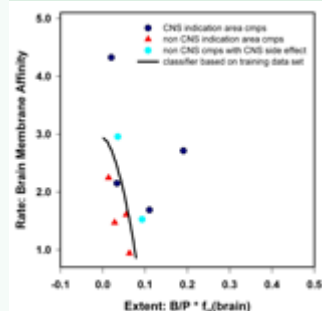
[ISBC](#) 2010, the 16<sup>th</sup> International Symposium on Bioluminescence and Chemiluminescence, is taking place in Lyon. The [NightSHADE](#) plant imaging system as well as the new Lumat<sup>3</sup> tube luminometer will be on display.

Berthold Technologies is happy to donate the [Marlene DeLuca prize](#) again to young scientists.

The NightSHADE plant imager will also be presented at the [EUCARPIA](#) Cereals Section of the National Institute of

Get in touch  
with Berthold  
Technologies

[www.  
Berthold.com  
/Bio](http://www.Berthold.com/Bio)



#### Upcoming Events:



Cambridge, UK  
April 6 - 8



Phoenix, AZ, USA  
April 12 - 15



Lyon, France  
April 19 - 23

[Find later events](#)

## Berthold on Facebook

Berthold Technologies has established a [Facebook page](#) with useful information about applications and products.

**Have a look and let us know your thoughts.**

## NightSHADE Plant Imager

Based on the experiences with the [NightOWL™](#) in vivo\* imager Berthold Technologies has developed the new [NightSHADE](#) imaging system with plant research in mind.

For applications such as gene research, circadian rhythms, regulation of plant growth, stress tolerance and drug screening the NightSHADE will offer unique features for plant research.

- ▶ LED-based plant illumination with daylight simulation
- ▶ 2<sup>nd</sup> camera for side-view imaging
- ▶ rotating table for 360° viewing

## New Application & Technical Notes

- ▶ Neutrophil Chemotaxis Assay performed on Twinkle LB 970

[Go to download page](#)

- ▶ Integration of Mithras and Centro plate readers with Hudson Robotics PlateCrane

[Go to download page](#)

[Have a look at other Application Notes](#)

## Recent Publications

- ▶ Measurement of Changes in Cytosolic Ca<sup>2+</sup> in Arabidopsis Guard Cells and Mesophyll Cells in Response to Blue Light; Plant and Cell Physiology 2009 50(2):360-373

[Go to abstract](#)

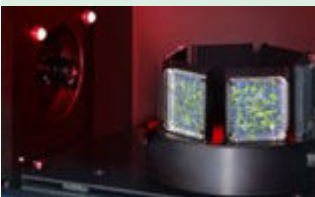
- ▶ Chimeric mouse tumor models reveal differences in pathway activation between ERBB family- and KRAS-dependent lung adenocarcinomas; Nature Biotechnology 28, 71-78 (2010)

[Go to abstract](#)

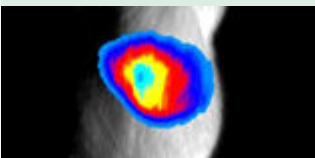
- ▶ Autophagy influences glomerular disease susceptibility and maintains podocyte homeostasis in aging mice; J Clin Invest. doi:10.1172/JCI39492

[Go to abstract](#)

[Have a look at all publications](#)



[Learn more about NightSHADE](#)



\* Some techniques for generating and/or detecting light in biological subjects are patented and may require licences from third parties. Users are advised to independently determine for themselves whether their activities infringe any valid patent.

[Unsubscribe from this newsletter](#)

**BERTHOLD TECHNOLOGIES GmbH & Co KG**

Headquarters/Germany +49 7081 177 200 • USA +1 865 483 1488 • UK +44 1582 76 1477

Japan +81 3 5825 3557 • Italy +39 03 9287 3064 • Switzerland +41 4 4871 2500

India +91 22 2765 5462 • BeNeLux +32 2251 6010 • France +33 1 3494 7900

China +86 10 8813 1974 • Austria +43 1 9142 2510 • Australia +61 3 9467 6277

Spain +34 91 361 4428 • Nordic +47 32 135 435 • Korea +82 31 776 0771

Please visit web page for other countries