

News September 2009



Get a TriStar with the purchase of a NightOWL™

60th Anniversary Promotion

60th Anniversary of Berthold

BERTHOLD TECHNOLOGIES is looking back to 60 years of company history. In the year 1949 Prof. Rudolf Berthold had laid the cornerstone for the company in the building of a former hotel. In the course of the past 60 years an enterprise with more than 300 employees and 3 business units has evolved. The innovative and high-quality measurement instruments are still developed and manufactured in Bad Wildbad, Germany, and sold via a worldwide network of affiliates and distributors.

Anniversary Promotions

BERTHOLD TECHNOLOGIES wants our valued customers to be able to participate in the celebrations. Until the end of 2009 a series of promotions will be issued each of them containing an attractive offer based around a specific instrument or a group of instruments. Details of all issued promotions can be found on our webpage www.Berthold.com/bio/promotions.

Stay tuned for the next newsletters and let your fellow researchers know about it!

And this is the present promotion:

Get a Tristar multimode reader with the purchase of a NightOWL™ in vivo imager

With the purchase of a NightOWL™ NC 100 in vivo imaging* system you will get a TriStar multimode plate reader in addition (see the applying terms and conditions below).

In many cases in vitro assays and cell culture experiments in microplates need to be performed to validate in vivo results. BERTHOLD TECHNOLOGIES is able to offer this unique combination of instruments – and in the course of the 60th anniversary of the company you will be able to benefit from a very attractive pricing by obtaining both instruments for the price of a NightOWL™ only.

[Check out the Anniversary Promotions](#)

Upcoming Shows:



**BIO
TECHNICA**
6. - 8. Oktober 2009
Hannover

We'll see you at:
Hall 9 Stand E02



**9th
IPNB
CONGRESS**
October 25-30
2009
St. Louis, Missouri, USA



MEDICA®
Weltforum der Medizin
Düsseldorf
18. - 21. Nov. 2009

[Learn more about NightOWL™](#)

Simply ask your local representative for a respective quotation.

NightOWL™ LB 983 in vivo imager

For in vivo imaging applications the NightOWL™ family is the solution meeting the demands of today's research:

- **Variable field of view due to moving camera with continuous height calibration**
- **High sensitivity due to deep cooled cameras with high quantum efficiency**
- **Many accessories for animal and plant research**
- **Luminescence and fluorescence applications**
- **Modern and intuitive Indigo™ software for acquisition and analysis**



[Learn more about TriStar](#)

The TriStar LB 941 multimode plate reader

is an attractive and compact instrument for microplate assays:

- **Luminescence/Fluorescence/Absorbance**
- **Dedicated Optical Path System (DOPS) for optimized conditions and best sensitivity**
- **JET injectors for most accurate and reliable reagent dispensing**
- **BRET and BRET² capability**



[Get in touch with Berthold Technologies](#)

www.berthold.com/Bio

Terms and Conditions

This promotional offer is valid on NightOWL™ LB 983 instruments.

The offer has no cash value.

The offer is valid with new instruments and new orders only.

The resulting overall discount must not exceed 25%.

No substitutions are allowed.

The offer is valid only for orders received by Berthold Technologies and its affiliated companies no later than December 4th 2009.

Berthold Technologies reserves the right to change the terms of or discontinue this promotion without further notice.

The offer is void where prohibited or restricted by law.

*** 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