

News October 2009



temperature control and
bottom reading for free

60th Anniversary Promotion

60th Anniversary of Berthold

BERTHOLD TECHNOLOGIES is looking back over 60 years of company history. In the year 1949 Prof. Rudolf Berthold had laid the foundation 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 supported via a worldwide network of affiliates and distributors.

Anniversary Promotions

BERTHOLD TECHNOLOGIES invites our valued customers 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 latest promotion:

Get the bottom reading module and the temperature control option for free with the purchase of a Twinkle fluorescence plate reader

This is a unique offer to extend the versatility of your plate fluorometer, e.g. for cell based assays.

Bottom reading can be an interesting instrument feature when measurements of adherent cells, with coloured buffer or with lids placed on top of the microplate, have to be performed to achieve the most efficient detection.

To have temperature control capabilities in an instrument is a must when enzymatic activities are monitored over time and when cellular assays are performed. Both a controlled and stable temperature during the whole

[Check out the Anniversary Promotions](#)

Upcoming Shows:



measurement time is important.

Get both at no additional cost
(see terms and conditions below).

Simply ask your local representative for
a respective quotation.

Twinkle LB 970 fluorescence plate reader

Is a dedicated fluorometer which has been designed
without compromise resulting in optimized performance
for fluorescent microplate assays.

- **low noise photomultiplier**
- **minimum scattering optics guarantee high sensitivity**
- **short and long term kinetics**
- **well scanning**
- **plate shaking**
- **dual excitation or emission**
- **bottom reading**
- **temperature control**
- **multiple plate formats incl. Petri dishes**

All these features guarantee the high sensitivity and
emphasise the versatility of the Twinkle fluorometer
for biochemical and cell based assays.

Terms and Conditions

This promotional offer is valid on Twinkle LB 970 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.

[Learn more
about
Twinkle](#)



[Get in touch
with Berthold
Technologies](#)

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

[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