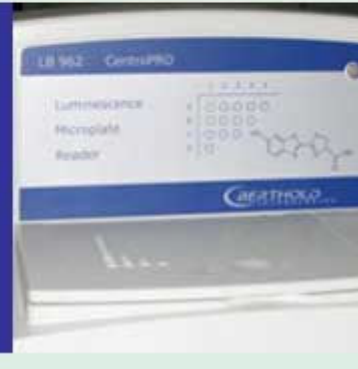


News August 2009



Win a Microplate Luminometer

60th Anniversary



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 headstone 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.

[Register for the Lottery to win a Microplate Luminometer](#)

Win a Microplate Luminometer

A lottery has been established within the anniversary activities to enable the valued customers of Berthold Technologies to participate in the celebrations. Until October 15th 2009 every interested scientist may simply register online via internet www.Berthold.com/bio/lottery.



After the registration will have been closed **3 microplate luminometers CentroPRO will be awarded**. On the occasion of **Biotechnica** exhibition 1 instrument will be awarded, too: registration may be done right at the Berthold Technologies **stand E02 (hall 9)**.

[Learn more about CentroPRO](#)

CentroPRO LB 962 Microplate Luminometer

CentroPRO is the newest member of the microplate luminometer family of Berthold Technologies. The instrument is characterized by

- ❑ **Easy operation via the innovative ICEware: wizard guided definition of parameters various export formats (Excel, PDF, text) intuitive menus**
- ❑ **High sensitivity due to Single Photon Counting technology and low noise detectors: less than 20 amol ATP less than 7.5 zmol Firefly luciferase**
- ❑ **Big dynamic range over 6 decades**



- ❑ **Small footprint with a width of only 30 cm**
- ❑ **All important glow-type luminescence applications**

CentroPRO is a universally applicable luminometer for research and routine use. It enables the measurement of all important glow-type luminescence applications: e.g. **reporter genes, dual reporter genes, ATP, cell viability, caspases, kinases, proteases, monoaminoxidase** or **Mycoplasma contamination**.

Terms and Conditions

This promotion is valid for CentroPRO LB 962.

The offer has no cash value.

No substitutions are allowed.

The promotion is valid until October 15th 2009.

The promotion is applicable to research organisations only.

The promotion is not applicable to Berthold Technologies employees and members of representing sales and service organisations.

The promotion is not applicable to employees and sales and service representatives of competing companies.

The promotion is not applicable for retail.

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.

The contact details collected will be distributed within and used by the Berthold Technologies sales organisations.



[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 1 9142 2510 • 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