

News September 2009



Make your instruments more versatile with injectors

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:

Pimp your instrument by adding JET injectors for reagent dispensing

Some applications like Calcium monitoring and enzyme kinetics require the addition of substrates or agonists immediately before or even during the measurement. Some labels show intrinsically fast reaction kinetics like Acridinium and Aequorin. Some detection chemistries are more sensitive in their flash versions, e.g. for ATP quantitation, and some detection chemistries are simply cheaper in their flash versions.

Many good reasons for having on-board reagent injectors.

Get the Berthold JET injectors free of charge and make

[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

your new instrument more versatile (see terms and conditions below).

Simply ask your local representative for a respective quotation.

JET injectors for reagent dispensing

The BERTHOLD TECHNOLOGIES JET injectors have a unique proprietary design based on Teflon bellows offering many advantages:

- **High accuracy (on average better than 99 % over entire volume range)**
- **High reproducibility (on average better than 99 % over entire volume range)**
- **Frictionless for longer lifetimes**
- **Cell-friendly**
- **Injection in reading position**

These instruments can be fitted with JET injectors:

- [Mithras LB 940](#) multimode plate reader
- [TriStar LB 941](#) multimode plate reader
- [Centro LB 960](#) plate luminometer
- [Centro XS³ LB 960](#) plate luminometer
- [Lumat LB 9507](#) single tube luminometer
- [Autolumat LB 953](#) multi tube luminometer

Terms and Conditions

This promotional offer is valid on Mithras LB 940, TriStar LB 941, Centro/Centro XS³ LB 960, Lumat LB9507 and Autolumat LB 953 instruments.

The offer is not valid for special configurations and OEM instruments.

The offer has no cash value.

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

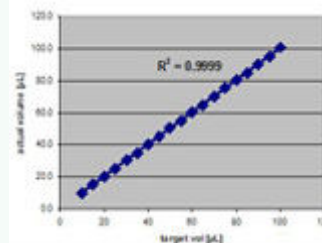
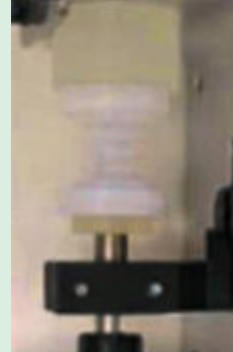
The resulting overall discount must not exceed 25%.

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 Microplate Readers](#)

[Learn more about Tube Luminometers](#)

[Get in touch with Berthold Technologies](#)

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