

News September 2009



Receive a set of cells with
the purchase of a FlowStar

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 corner stone 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:

Receive up to 2 measuring cells free of charge with the purchase of a FlowStar radioactivity monitor for HPLC systems

A unique offer to configure your ideal radio HPLC system.

With the purchase of a FlowStar radioactivity monitor you can choose between a variety of cells and accessories to be added free of charge to complete your dedicated HPLC system (see terms and conditions below):

- set of special UPLC cells
- set of regular solid cells
- set of gamma cells
- set of admixture cells for 3H applications

[Check out the Anniversary Promotions](#)

Upcoming Shows:



Simply ask your local representative for a respective quotation.

FlowStar LB 513 radio HPLC monitor

Radiolabelled compounds for drug, pesticide, metabolism and pharmacokinetic studies provide very high sensitivity and specificity which is not available with mass detection, UV, fluorescence or conductivity measurements.

The FlowStar represents the latest enhancement in flow-through detectors for Radio HPLC:

- **Highly sensitive detection unit operated in coincidence counting mode**
- **Automatic detection and identification of measuring cells**
- **Touch screen for easy operation**
- **Multiple input and output connectors**
- **RadioStar PC software**

The system is completed by the application specific measuring cells:

- Liquid Scintillation (admixture) cells
- Solid Scintillator cells
- Microbore cells
- LC/MS cells
- Fast LSC cells (UPLC, UHPLC, RLC)
- Preparative SFC cells
- Cerenkov cells for high energy beta emitters
- Gamma Isotopes and Positron Emitter (PET) cells
- Special Iodine cells
- GLP/GMP validation test cells
- Customized cells

Terms and Conditions

This promotional offer is valid on FlowStar LB 513 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 FlowStar](#)



[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