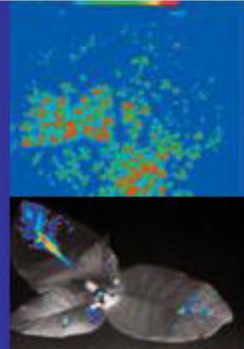




News October 2009

## NightSHADE

A new concept for  
plant research



### Newsletter Topics

- ❑ IPMB 2009 in St. Louis, MO, Oct. 25<sup>th</sup> - 30<sup>th</sup>
- ❑ NightSHADE - the new dedicated *in vivo* plant imager
- ❑ 60<sup>th</sup> anniversary promotions of Berthold Technologies

### 60<sup>th</sup> Anniversary of Berthold

Berthold Technologies is looking back over 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.

Until the end of 2009 a series of promotions will be issued. Details of all issued promotions can be found on our webpage [www.Berthold.com/bio/promotions](http://www.Berthold.com/bio/promotions).

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

### NightSHADE dedicated *in vivo* plant imager

Based on our experience with the NightOWL *in vivo*\* imager, Berthold Technologies has developed the new NightSHADE imaging system with plant research in mind.

For applications such as gene research, circadian rhythms, regulation of plant growth, stress tolerance and drug screening the NightSHADE will offer unique features for plant research.

- ❑ Top-view emCCD camera with high resolution
- ❑ LED-based plant illumination with daylight simulation
- ❑ 2<sup>nd</sup> camera for side-view imaging
- ❑ x-y table and rotating table for 360° viewing
- ❑ temperature & humidity control



[Check out the Anniversary Promotions](#)

Upcoming Shows:



**MEDICA®**

Weltforum der Medizin

Düsseldorf  
18. - 21. Nov. 2009

[Learn more about NightSHADE](#)

**IPMB 2009, St. Louis, MO, Oct. 25<sup>th</sup> - 30<sup>th</sup>**

Visit the exhibition and have a look at Berthold Technologies solutions for plant research:

**NightSHADE**

in vivo imager



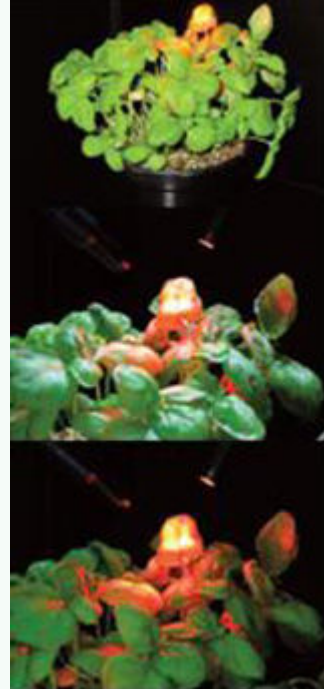
**Centro**

microplate reader



**Luminescence**

flow through



**Terms and Conditions**

This promotional offer is valid on instruments as indicated on the webpage.

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.**

[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