

Newsletter September 2010

in vivo Plant Imaging issue

Newsletter Topics

- ▶ **New Application Note: GFP**
- ▶ **Berthold Technologies at WMI, ILMAC and Cancer Association Meeting**
- ▶ **Recent Publications**

New Application Note

In vivo imaging of plants became a common research tool all over the world for gene expression, studies of circadian rhythm, etc. Plants are transfected either with Luciferase or photoproteins like GFP, YFP or dsRED.

In this application note we could show that excitation of GFP-transfected plants in Petri dishes filled with agar can be easily performed by trans-illumination, without reflection on the plastic material as is usually the case with epi-illumination.

- ▶ Trans-illumination – a Solution for Excitation of GFP Transfected Plants in Petri Dishes for *In Vivo* Imaging

[Go to download page](#)

[Have a look at other Application Notes](#)

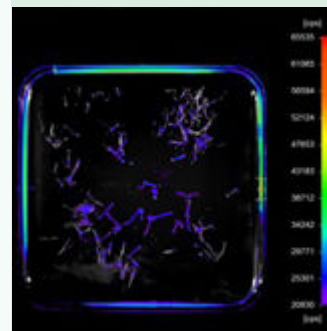
Berthold Technologies present at WMIC, ILMAC and Cancer Association Meeting

The **World Molecular Imaging Congress 2010** takes place in Kyoto, Japan, September 8-11th. Have a look to the imaging solutions from Berthold Technologies. We will welcome you on Stand No. 423!

The **ILMAC 2010**, Industry Exhibition for Research & Development in Pharmaceuticals, Chemicals and Biotechnology will be held September 21-24th in Basel, Switzerland. Berthold Technologies will present its **new Lumat³ tube luminometer**. Please come to Hall 1.1 - Stand A62!

Get in touch
with [Berthold Technologies](#)

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



GFP-transfected *Arabidopsis thaliana* plants trans-illuminated with 315 nm and measured in NightSHADE.



[Learn more about NightSHADE](#)

Upcoming Events:



Kyoto, Japan
September 8 - 11

ILMAC
21. - 24.09.2010

Basel, Switzerland

Visit us at the **69th Annual Meeting of the Japanese Cancer Association** in Osaka, Japan, September 22-24th.



Osaka, Japan
September 22 - 24

[Find later Events](#)

Recent Publications

- ▶ Protein Kinase C-induced Phosphorylation of Orai1 Regulates the Intracellular Ca⁺⁺ Level via the Store-operated Ca⁺⁺ Channel; The Journal of Biological Chemistry, 285, 25720-25730

[Go to abstract](#)



Calcium Assay with Fura-2 measured with Mithras LB 940

- ▶ Cell turnover in tissues of the long-lived ocean quahog *Arctica islandica* and the short-lived scallop *Aequipecten opercularis*; Marine Biology Volume 157, Number 6, 1283-1292

[Go to abstract](#)



Caspase Assay was measured in TriStar LB 941 - samples from various clam species!

- ▶ Effects of the Combination of Favipiravir (T-705) and Oseltamivir on Influenza A Virus Infections in Mice; Antimicrobial Agents and Chemotherapy 2010 January; 54(1): 126-133.

[Go to abstract](#)

[Have a look at all publications](#)



Viral neuraminidase inhibition assay (NA-Star detection kit, Applied Biosystems) was measured with Centro LB 960

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.

[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