

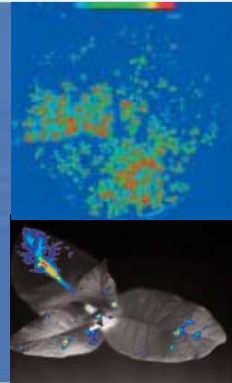
detect and identify



NEWS June 2009

NightSHADE

a new concept for Plant Research



Newsletter Topics

- Berthold Technologies at AgriGenomics 2009 in London
- NightSHADE – the new dedicated *in vivo* plant imager

NightSHADE – the dedicated *in vivo* plant imager

Based on the experiences 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
- 2nd camera for side-view imaging
- x-y table and rotating table for 360° viewing
- temperature & humidity control

AgriGenomics 2009, London, 2nd/3rd July



- Attend the [Workshop in vivo plant imaging](#).
- Listen to the *Technology Spotlight* [NightOWL in vivo imaging system](#) covering plate reader applications as well
- Visit the exhibition to see Berthold Technologies solutions for plant research:

NightSHADE
in vivo imager



Centro
microplate reader



Luminescence
flow through



[Learn more
about the
NightOWL
in vivo
imager](#)



[Get in touch
with Berthold
Technologies](#)

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

[Unsubscribe from this newsletter](#)