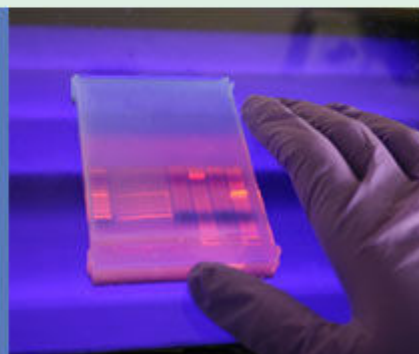


NEWS July 2009

New Imaging Systems for Gels & Blots



Newsletter Topics

- ❑ **Berthold Technologies has started to distribute the approved DNR gel and blot imaging systems**
- ❑ **Introduction of the new MiniLumi Gel Doc system**
- ❑ **The complete range of imaging system for gels and blots**

DNR imaging systems available through Berthold Technologies

In January 2009 Berthold Technologies has signed an agreement with DNR Bio-Imaging Systems to distribute DNR's imaging solutions for gels and blots stained with coloured, fluorescent or luminescent markers. DNR has a long tradition of developing and manufacturing scientific imaging systems. All systems are characterized by sophisticated illumination and camera types.

Berthold Technologies Germany and its offices in BeNeLux, France, Italy, Switzerland, UK, USA, India and Japan are now your qualified and reliable partners for sales and support of DNR imaging systems.

MiniLumi

The MiniLumi is a brand new system for fluorescent and colorimetric gel documentation with 1-click image capturing. Typical applications include Ethidium bromide, Coomassie® Brilliant Blue, SYPRO™ Orange, SYPRO™ Ruby, Silver stain and Q-Dots. With outstanding features like

- ❑ 2 Megapixel CCD camera
- ❑ 16 bit TIF files
- ❑ F/1.2 lens
- ❑ Trans-UV and trans-whitelight illumination
- ❑ Epi-UV and epi-whitelight illumination
- ❑ Firewire interface

the MiniLumi offers high quality at affordable cost.

[Learn more about DNR imaging systems](#)

[Download application notes about gel imaging](#)



MiniBIS Pro

The state-of-the-art and extremely flexible gel documentation system is characterized by a super-bright lens, a dual transilluminator with interchangeable drawers and the possibility to take the camera part off the system to put it onto separate transilluminators.

MiniBIS Pro features are:

- ❑ 1.3 Megapixel CMOS camera
- ❑ 16 bit TIF files
- ❑ F/0.95 lens
- ❑ Field of view up to 21 cm
- ❑ Trans-UV and trans-whitelight illumination

LumiBIS

The LumiBIS introduces full automation into gel documentation by its smart dark chamber with motorized camera movement to change the field of view, motorized zoom and focus. In addition to various standard trans-illumination and epi-illumination options unique multi-LED illumination may be used extending the range of applications by SYBR™ Green, ProQ™ Diamond, eGFP, eYFP, dsRed, Alexa594 and many more. LumiBIS features include:

- ❑ 2 Megapixel CCD camera
- ❑ 16 bit TIF files
- ❑ F/1.2 zoom lens, 8 - 48 mm
- ❑ Trans-UV and trans-whitelight illumination
- ❑ Epi-UV and epi-whitelight illumination
- ❑ LED illumination: 470, 500, 525, 568, 588, 621 nm

MF ChemiBIS

The superior system for gel and blot imaging and analysis is equipped with a high resolution low noise cooled CCD camera enabling the detection of luminescent Western blots, e.g. ECL™, Super Signal™, CDPStar™ or CSPD™. Highest degree of automation through motorized camera, zoom and focus and multiple illumination sources underline the flexible usage for an almost unlimited range of applications with colorimetric, densitometric, fluorescent and luminescent detection. NLD TL-100 software is used for image analysis. MF ChemiBIS features:

- ❑ 3.2 Megapixel CCD camera
- ❑ Deep cooling system (-70 °C)
- ❑ High full well capacity
- ❑ F/0.95 super bright lens
- ❑ 16 bit TIF files
- ❑ Trans-UV and trans-whitelight illumination
- ❑ Epi-UV and epi-whitelight illumination
- ❑ LED illumination: 470, 500, 525, 568, 588, 621 nm



[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 1 9142 2510 • France +33 1 3494 7900