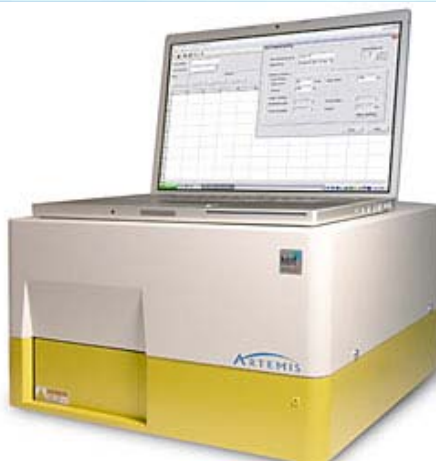


Artemis

HTRF[®] Microplate Reader



Berthold Technologies completes HTRF[®] reader portfolio

Starting August 15th, 2008, BERTHOLD TECHNOLOGIES, a leading manufacturer of microplate readers for diagnostics, research and drug discovery has taken over the worldwide distribution of Furuno's Artemis HTRF[®] microplate reader.

The Artemis K-101 has been designed for optimum performance with CISbio's HTRF[®] reagents. The HTRF[®] approved instrument is extremely easy to use and guarantees superior performance even with demanding applications.

- **High energy excitation**
- **Low background optics**
- **Dedicated filters**
- **Wide dynamic range**
- **High sensitivity**
- **HTRF[®] ratio calculation**

The Artemis completes the portfolio as Berthold Technologies is now able to offer the most extensive range of microplate readers with TRF and HTRF[®] functionality

| Instrument | Reading Technologies |
|---------------------------|---|
| Artemis | HTRF [®] |
| Mithras HTRF Basic | Luminesc., Fluoresc., TRF, HTRF [®] |
| Mithras HTS II | Luminesc., BRET, Fluoresc., VIS Abs., Fluor. Polar., AlphaScreen™, TRF, HTRF [®] |
| Mithras HTS III | Luminesc., BRET, Fluoresc., UV/VIS Abs., Fluor. Polar., TRF, HTRF [®] |

[Learn more about Artemis](#)

[Learn more about Mithras](#)



htrf

[Get in touch with Berthold Technologies](#)

www.Berthold.com/Bio