

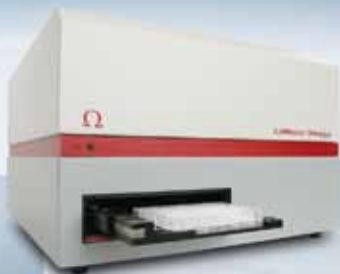
LUMIstar Omega



The Dedicated Luminescence Microplate Reader

- High performance microplate reader for luminescence
- Measurement at up to eight distinctive wavelengths
- Top and bottom reading
- BRET, optional with Simultaneous Dual Emission detection
- Up to two reagent injectors
- Upgradeable to FLUOstar Omega or POLARstar Omega including multiple detection modes and UV/Vis spectrometer
- Powerful MARS Data Analysis Software





LUMIstar Omega - Technical Specifications

Due to the modularity of BMG LABTECH's instruments, all or combinations of the features below can be installed at purchase or upgraded at any time. Please contact your local representative for more details or a quote.

Detection Modes	Luminescence (flash and glow) - including BRET
Measurement Modes	Top and bottom reading Endpoint and Kinetic measurements Sequential Multi Emission measurements Simultaneous Dual Emission measurements Ratiometric measurements
Microplate Formats	6 to 384-well plates, user-definable
Detectors	Side window, current type photomultiplier tube
Optical Filters	One emission filter wheel for 8 filters
Spectral Range	240 - 740 nm
Sensitivity	20 amol/well ATP DLR certified
Read Times	Flying mode: 20 s (96), 55 s (384)
Reagent Injections	Up to 2 built-in reagent injectors Injection at measurement position (6 to 384-well) Individual injection volumes for each well (3 to 350 µL) Variable injection speed up to 420 µL / s Up to four injection events per well Reagent back flushing
Shaking	Linear, orbital, and double-orbital with user-definable time and speed
Gas Vent	System to inject an atmosphere or to pull a vacuum into the reader
Incubation	+5°C above ambient up to 45°C or 60°C
Software	License-free software package including Reader Control and MARS Data Analysis Software
Dimensions	Width: 44 cm, depth: 48 cm, height: 30 cm; weight: 28 kg
Accessories	
Stacker	Magazines for up to 50 plates - continuous loading feature
THERMOstar	Microplate Incubator and Shaker
Filters	Optimized for dyes, fluorophores and specific assays Filters for all applications from UV to NIR Customized filters available upon request
Upgrades	Upgrades to include options such as additional detection modes, reagent injectors, extended temperature control, etc. are available. Please contact your local representative for more information.



Headquarters

Germany
BMG LABTECH GmbH
Allmendgrün 8
77799 Ortenberg
Tel. +49 781 96968 -0
Fax +49 781 96968 67
germany@bmglabtech.com

Australia

BMG LABTECH Pty. Ltd.
2/24 Carbine Way
Mornington, Victoria, 3931
Tel. +61 3 5973 4744
Fax +61 3 5973 4711
australia@bmglabtech.com

France

BMG LABTECH SARL
7, Rue Roland Martin
94500 Champigny s/Marne
Tel. +33 1 48 86 20 20
Fax +33 1 48 86 47 07
france@bmglabtech.com

Japan

BMG LABTECH JAPAN Ltd.
2F TS-1 Building
1-6-2, Shimo-cho
Omiya-ku
330-0844 Saitama City
Tel. +81 48 647 7217
Fax +81 48 647 7218
japan@bmglabtech.com

UK

BMG LABTECH Ltd.
5 Merlin Centre
Gatehouse Close
Aylesbury
HP19 8DP
Tel. +44 1296 336650
Fax +44 1296 336651
uksales@bmglabtech.com

USA

BMG LABTECH Inc.
13000 Weston Parkway
Suite 109
Cary, NC 27513
Tel. +1 877 264 52 27
Fax +1 919 678 16 40
usa@bmglabtech.com