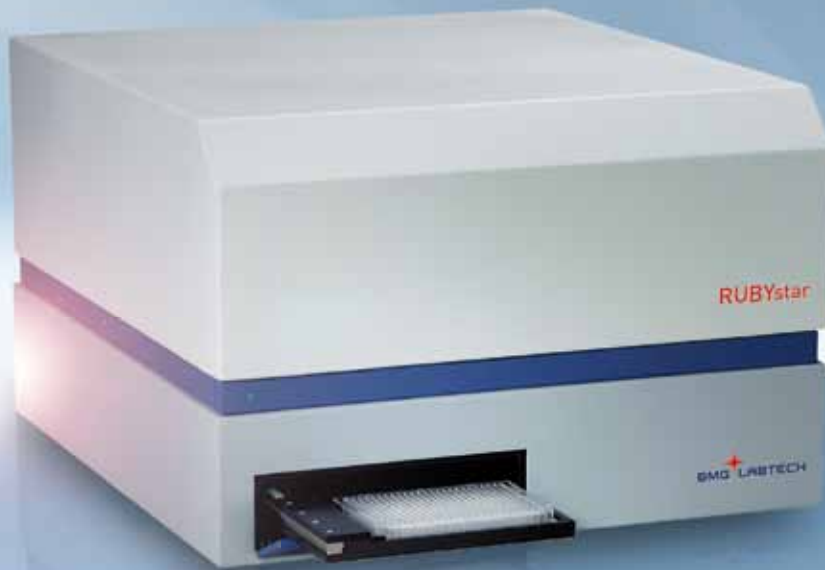


RUBYstar



High-performance Laser-based HTRF[®] Microplate Reader



- Time-resolved fluorescence and HTRF[®]
- High energy nitrogen laser
- Single photon counting
- Fastest read times with Simultaneous Dual Emission detection
- Unique decay curve monitoring with 2 μ s resolution

www.bmglabtech.com

**BMG LABTECH**

The Microplate Reader Company

RUBYstar - 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	Homogeneous Time-Resolved Fluorescence
Measurement Modes	Top Reading Ratiometric measurement with real-time ratio
Microplate Formats	Up to 384-well plates
Light Source	High-energy nitrogen laser (337 nm)
Detector	Lens based dual channel photomultiplier system for Simultaneous Dual Emission at 620 nm and 665 nm Single photon counting
Read Times	5 flashes 55 s (96), 170 s (384) 20 flashes 130 s (96), 475 s (384)
Software	License-free Software package including Reader Control and Data Evaluation Software
Dimensions	Width: 44 cm, depth: 48 cm, height: 30 cm; weight: 29 kg
Accessories	
Stacker	Magazines for up to 50 plates - continuous loading feature
THERMOstar	Microplate Incubator and Shaker
Upgrades	Upgrades to include options such as reagent injectors 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
Morningside, 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

HTRF is a registered trademark of Cisbio International.

Limit of detection was calculated according to the IUPAC standard: $3(SD_{blank})/slope$
Specifications are subject to change without notice.

www.bmglabtech.com