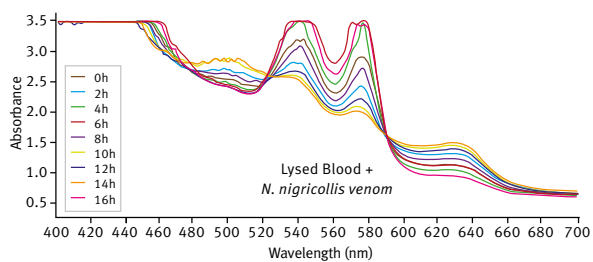




Microplate reader with ultrafast UV/Vis spectrometer

The SPECTROstar® Omega is a dedicated absorbance microplate reader that can capture either a full-absorbance spectrum or up to eight discrete wavelengths simultaneously. Upgrades can transform the SPECTROstar Omega into a powerful multi-mode reader with up to six detection modes, Simultaneous Dual Emission and injectors.

- Captures UV/Vis absorbance spectra from 220-1000 nm
- Spectrum acquisition faster than 1 second per well
- Plate formats up to 1536-well
- Plate shaking and incubation up to 45°C or 65°C
- Up to two integrated reagent injectors
- Upgradeable to powerful multi-mode reader
- Compatible with LVis plate and Stacker



Time-dependent change in haemoglobin absorbance spectrum in the presence of *N.nigricollis* venom. Lysed sheep blood incubated with PBS only or *N. nigricollis* venom. Spectrum recorded at 2 hour intervals.

SPECTROstar® 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	UV/Vis absorbance	
Measurement modes	Endpoint and kinetic measurements Well scanning	
Microplate formats	6 to 1536-well plates, user-definable	
Light source	High energy xenon flash lamp	
Detectors	Spectrometer with CCD	
Spectral range	Absorbance spectrometer: 220 to 1000 nm	
Sensitivity	ABS with spectrometer	Spectral range: 220 to 1000 nm Full spectrum captured in < 1 s/well Selectable spectral resolution: 1 - 10 nm OD range: 0 to 4 OD Accuracy: < 1% at 2 OD Precision: < 0.5% at 1 OD and < 0.8% at 2 OD Reproducibility: ±0.005 OD for 0 - 2 OD range
Read times	Dependent on test definition	
Reagent injection	Up to 2 built-in reagent injectors Injection at measurement position (6 to 384-well) Individual injection volumes for each well (3 to 500 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	+4°C above ambient up to 45°C or 65°C	
Software	Multi-user Reader Control and MARS Data Analysis Software package included	
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	
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.	

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

© 2017 All rights reserved. All logos and trademarks are the property of BMG LABTECH.

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

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

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

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

UK
BMG LABTECH Ltd.
5 Alton House Office Park
Gatehouse Way
Aylesbury
HP19 8XU
Tel. +44 1296 336650
uksales@bmglabtech.com

USA
BMG LABTECH Inc.
13000 Weston Parkway
Suite 109
Cary, NC 27513
Tel. +1 877 264 5227
usa@bmglabtech.com

www.bmglabtech.com