

# Evaluation Plate



## Functional validation of your plate reader

The Evaluation Plate verifies and validates the optical functionality of all BMG LABTECH plate readers for quality control and/or qualification purposes. Simplified fluorescence intensity, luminescence and absorbance qualification tests save time and avoid the use of liquid chemistry.

The Evaluation Plate is equipped with:

- Reference optical filters
- Solid fluorescence standards
- LEDs with defined emission intensities

Test measurements, data reductions and attainment/failure reports are automatically performed by the software.

## Tests and compatible instruments

Here below you can find the qualification features divided by reader

Mode	Parameter	CLARIOstar	Omega	PHERAstar
Absorbance	Absolute accuracy	●	●	●
	Wavelength accuracy	●	●	●
	Reproducibility	●	●	●
	Linearity	●	●	●
Fluorescence intensity	Linearity	●	●	●
	Reproducibility	●	●	●
	Wavelength scan	●		
Luminescence	Linearity	●	●	●
	Reproducibility	●	●	●
	Intensity	●	●	●
	Wavelength scan	●		

### Headquarters Germany

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

### Australia

BMG LABTECH Pty. Ltd.  
2/24 Carbine Way  
Morningside, 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



Made in Germany