

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.
8 Bell Business Park
Smeaton Close
Aylesbury
Bucks
HP19 8JR
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