



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 standards.

The Evaluation Plate is equipped with:

- NIST-traceable absorbance reference filters
- Solid fluorescence standards
- Solid luminescence standards

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

Validation tests and compatible readers:

Mode	Absorbance				Fluorescence intensity			Luminescence			
Parameter	Absolute accuracy	Wavelength accuracy	Reproducibility	Linearity	Linearity	Reproducibility	Wavelength scan	Linearity	Reproducibility	Intensity	Wavelength scan
CLARIOstar <i>Plus</i>	●	●	●	●	●	●	●	●	●	●	●
Omega	●	●	●	●	●	●		●	●	●	
PHERASTAR <i>FSX</i>	●	●	●	●	●	●		●	●	●	
SPECTROstar <i>Nano</i>	●	●	●	●							

For further information on compatibility and technical requirements, please contact your local sales representative.

www.bmglabtech.com