

LVis Plate



Measure low-volume samples with ease!

BMG LABTECH's proprietary LVis Plate is a low-volume microplate that incorporates a cuvette slot and optional performance testing features. The LVis Plate is ideal for quick and easy low-volume absorbance measurements of DNA, RNA, protein samples, or spectral scanning.

The LVis Plate has the following outstanding features:

- Sixteen microdrop well sites for 2 μ L samples
- A horizontal position for a standard cuvette
- Four NIST-traceable optical density filters for quality control assessment
- A holmium oxide filter to assess wavelength accuracy



Easy pipetting of up to sixteen 2 μ L samples using the integrated multichannel pipette tip support.

The microdrop well sites are easily accessible. Left- and right-handed tip supports for 8 channel pipettes make pipetting multiple samples simple and easy.

The quick-clean surface easily wipes clean to prepare for further measurements.

The LVis Plate is compatible with all UV/Vis absorbance spectrometer-based BMG LABTECH microplate readers.

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
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



Made in Germany