

DATA SHEET

SpectraTest ABS2 Validation Plate

Comprehensive, industry-standard absorbance microplate reader validation



The SpectraTest® ABS2 Validation Plate is our next generation, automated, absorbance validation tool for Molecular Devices microplate readers.

Extended linear dynamic range

Test up to 3.0 OD for our latest microplate readers, SpectraMax® iD3, iD5, ABS, and ABS Plus readers, and up to 2.0 OD for earlier models. Validating the absorbance optical performance of your microplate reader is as simple as reading a plate.

Expanded wavelength coverage

Validate the optical performance of your absorbance reader in the visible range between 405–750 nm with NIST and NMI traceability. UV validation services are available upon request.

Automated certificate value entry

Reduce potential errors in the process of entering certificate values using our quick and easy copy/paste process. Values from the digital certificate are copied to the validation protocol and verified with a checksum to further increase process security.

Accredited recertification service

Recertify your ABS2 plate at our ISO 17025 accredited* facility in San Jose, California. Plates are handled by certified validation plate specialists and new certificate values for your plate are produced.

*View our [ISO 17025 certificate of accreditation](#).

KEY FEATURES

- Auto-entry for certificate values
- Preconfigured validation protocols
- NIST and NMI traceability
- Recertification services
- StakMax® Microplate Handling System compatible

Tests performed with the ABS2 plate

The ABS2 plate validates your reader using the following tests:

- Endpoint and kinetic baseline noise
- Optical alignment
- Ultimate dark
- Photometric accuracy (linearity)
- Photometric precision (reproducibility)
- Stray light
- Wavelength accuracy
- Wavelength precision

Description	Details	Part number
SpectraTest ABS2 Absorbance Validation Plate (Requires SoftMax® Pro Software version 6.5.1 or higher.)	Compatible with SpectraMax iD3, iD5, i3x, i3, M2, M2e, M3, M4, M5, M5e, Plus 384, ABS, ABS Plus, FlexStation 3 readers	0200-6191
Gray replacement validation plate case	Compatible with SpectraTest ABS2, FL1, LM1 plates	5074718
SpectraTest recertification		Part number
SpectraTest ABS2 Plate re-cert		S9200-0186

Contact Us

Phone: [+1.800.635.5577](tel:+18006355577)
 Web: www.moleculardevices.com
 Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada	+1.800.635.5577	China (Beijing)	+86.10.6410.8669	Japan	+81.3.6362.9109
United Kingdom	+44.118.944.8000	China (Shanghai)	+86.21.3372.1088	South Korea	+82.2.3471.9531
Europe*	00800.665.32860	Hong Kong	+852.3971.3530		

*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland