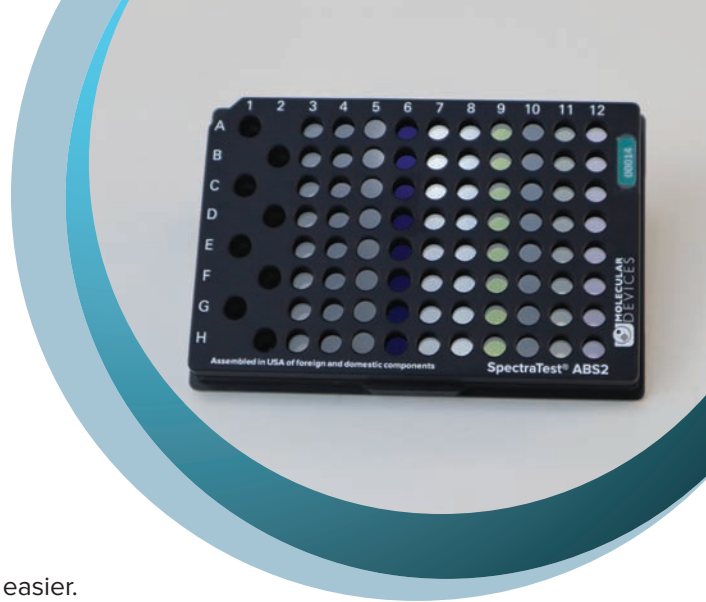


Upgrade to the next generation SpectraTest ABS2 Absorbance Validation Plate



Validating the optical performance of your plate reader has never been easier. The new SpectraTest® ABS2 Absorbance Validation Plate simplifies the certificate entry process to reduce human error and expands the testing capabilities to include more wavelengths and an extended linear dynamic range. Enhanced test methods and vendor recommended acceptance criteria ensure your reader is thoroughly tested against specifications we consider critical to achieving quality results.

Feature	ABS1 Validation Plate	ABS2 Validation Plate
Linear dynamic range (photometric accuracy and precision tests)	Up to 1.7 OD: all supported readers	Up to 3.0 OD: SpectraMax® iD3, iD5, ABS, ABS Plus readers Up to 2.0 OD: SpectraMax i3/i3x, M Series, Plus 384, and FlexStation® 3 readers
Wavelengths tested for photometric accuracy and precision (NIST and NMI traceable)	440, 465, 546, 590, 635 nm	405, 440, 465, 490, 546, 590, 635, 650, 750 nm
Wavelength accuracy test range	330*–680 nm	330*–775 nm
Certificate of calibration	Paper-based	Digital and paper-based
ISO 17025 accredited recertification service	•	•
8+ tests to validate optical performance	•	•
8 replicates of each filter type tested	•	•
Simple to understand test calculations and results		•
Improved test methods and statistical significance of results		•
Overall acceptable/out of specification result		•
Automated EZinCert certificate value entry		•
Automation friendly design for use with StakMax® Microplate Handling System		•
Updated packaging for easier shipping and handling		•
ABS2 validation plate		Part number
ABS2 validation plate (requires SoftMax® Pro Software version 6.5.1 or higher)		0200-6191
ABS2 recertification service		S9200-0186
Gray replacement case		5074718

*Validated wavelengths are dependent on the reader model.

Contact Us

Phone: +1.800.635.5577
 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