

SpectraDrop Micro-Volume Microplate

A micro-volume solution for benchtop microplate readers

KEY FEATURES

- **Elegant simplicity with multi-user design**
- **User versatility with different volume size and density options**
- **Cleaning convenience with hassle-free handling**
- **Unmatched throughput with up to 64 samples per plate**

Formerly known as the μ Max, Molecular Devices' unique SpectraDrop™ Micro-Volume Microplate offers the highest throughput solution for low volume measurement available on the market today. Innovative and flexible design features enable accelerated sample preparation time and increased laboratory productivity of DNA, RNA, and protein samples as low as 2 μ L. The SpectraDrop Micro-Volume Microplate assures uniform and reproducible analysis and integrates seamlessly with the StakMax® Microplate Stacker for greater research capacity.

Elegant simplicity

The SpectraDrop Microplate incorporates a specially designed adapter and a slide pair providing a uniform multi-sample assembly. The simplicity of this arrangement allows each user to have their own slide combination, making it ideal for multiuser laboratories.

The SpectraDrop Microplate Adapter has two side pockets for easy slide access and tight slide tolerances for a no-smear experience. The bottom slide is Teflon-

coated to ensure precise droplet placement, while the top slide has side spacers for 0.5mm (2 μ L) or 1mm (4 μ L) pathlength and evaporation control, making it a uniform and reproducible data analysis tool.

User versatility

The SpectraDrop Microplate comes in two density configurations, 24-wells per plate and 64-wells per plate, and two volume options, 2 or 4 μ L, allowing users to select the configuration ideal for their needs.

Cleaning convenience

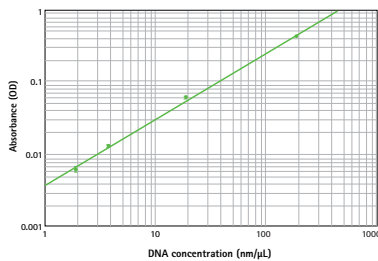
The two-slide design makes the cleaning process a breeze. The users can quickly wipe each slide, autoclave them, sonicate or simply replace them without having to worry about any calibration steps.

Unmatched throughput

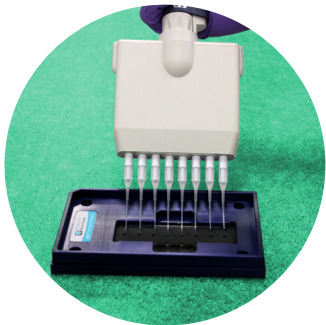
With up to 64 wells per sample and compatibility with Molecular Devices StakMax Microplate Stacker, the SpectraDrop Micro-Volume Microplate offers the highest low volume throughput available on the market today.



Unique design. The SpectraDrop Micro-volume microplate consists of specially designed adapter, Teflon coated bottom slide, and top slide with evaporation reducing spacers.



Reliable results. SpectraDrop Microplates are supported by SoftMax® Pro Microplate Data Acquisition and Analysis Software, which offers ready-to-read protocols.



Simplicity of use. For simplicity of use, the slide design is compatible with 8- and/or 16-channel pipettors.

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

Technical specifications

Sensitivity	2ng/μL (dsDNA)*
Uniformity	< 5% CV
24-well slide	Compatible with 8-channel pipettors
64-well slide	Compatible with 8- or 16-channel pipettors

Ordering information

Part number	Description	Details
0200-6262	SpectraDrop Starter Kit	Contains one microplate adapter, two 24-well low volume bottom slides, and one of each cover slide (2μL and 4μL)
0200-6267	SpectraDrop High-Throughput Screening (HTS) Kit	Contains one microplate adapter, five 24-well low volume slides, five 64-well low volume bottom slides, and five of each cover slide (2μL and 4μL)
0200-6263	SpectraDrop Micro-Volume Refills, 24-Well Bottom Slides	Three low volume 24-well bottom slides
0200-6264	SpectraDrop Micro-Volume Refills, 64-Well Bottom Slides	Three low volume 64-well bottom slides
0200-6265	SpectraDrop Micro-Volume Refills, 2μL Cover Slides	Three 2μL top slides (0.5 mm clear path length)
0200-6266	SpectraDrop Micro-Volume Refills, 4μL Cover Slides	Three 4μL top slides (1.0 mm clear path length)

* dsDNA sensitivity specified for the following instruments: SpectraMax® 190 Readers, SpectraMax® Plus384 Readers, SpectraMax® M-Series Readers, SpectraMax® Paradigm® Systems, FlexStation® 3 Systems