

## StakMax Microplate Stacker

Walk-away automation for benchtop microplate readers

### Key Features

- Integrates with SpectraMax® Readers
- Robust and reliable
- Easy to set up and use
- Flexible operation
- Leverages the power of SoftMax Pro Software

A key factor in achieving higher throughput in laboratory experiments is the ability to easily process batches of microtiter plates. Options for automating this process are often expensive, hard to use, or problematic. To address these roadblocks, Molecular Devices® has developed the StakMax® Microplate Stacker. The StakMax System is an easy-to-set-up, powerful, and reliable stacking solution for microplate readers from Molecular Devices. The system is operated through the industry-leading SoftMax® Pro Microplate Data Acquisition and Analysis Software and provides an easy-to-use interface for both novice and advanced users alike. The StakMax Stacker provides options for 20-, 40-, and, 50-plate magazines, and barcode reading.

### Simple to use

The integration of the StakMax System with a microplate reader is simple and user-friendly. All the tools necessary for setup come standard with the system, including a software calibration procedure. The entire setup process takes approximately 15 minutes and, once the system is aligned, it is ready to operate. The StakMax Stacker is programmed through SoftMax Pro Software for easy setup of batches or single-plate operation.

### Flexible operation

The StakMax Stacker offers flexibility in configuring and operating the system. The system comes with a single plate loading adapter and a choice of magazines for 20, 40 or 50 SBS-standard 96- and 384-well microplates. Users can choose to process all plates within the magazine or select a specific number. The StakMax Stacker also allows users to feed plates into the loading area manually.

A barcode reader accessory can be purchased with the system or added later and is easily installed onsite. The barcode reader can be configured to read either the short or long side of the microplate.

### Leverages SoftMax Pro Software

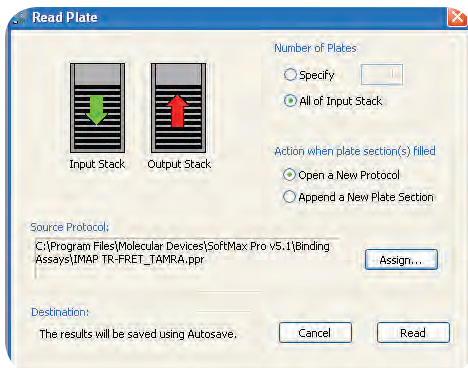
All of the StakMax system's configuration options are controlled through the standard SoftMax Pro Data Analysis Software. The system is also compatible with SoftMax Pro GxP Software, enabling bioassay validation in a secure FDA 21 CFR Part 11, GLP-/GMP-compliant environment.

## Flexible integration options



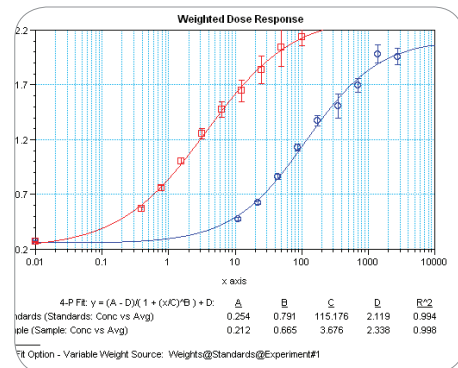
Integrate SpectraMax Absorbance, Fluorescence, Luminescence, and Multi-Mode Readers. System setup and calibration are controlled from within SoftMax<sup>®</sup> Pro Software.

## Easy batch setup



StakMax System Control Window: With a few mouse clicks, users can set up a batch process, all using existing SoftMax Pro Software protocols.

## Powerful data analysis tools



Complete data analysis for multi-plate assays and over 120 assay protocols are included with SoftMax Pro Software. Optional GxP compliance tools available.

## Reliable, robust automation

Molecular Devices has over 25 years of experience working with automation in the life sciences industry. We have leveraged this knowledge to develop a robust, reliable solution for automating microplates. With its superior reliability, users can set up the StakMax System for "walk-away" batch processing, which means less hands-on time processing plates.

## Compatibility

The StakMax System requires SoftMax Pro Software version 5.1 or greater and is compatible with the following benchtop microplate readers:

- SpectraMax<sup>®</sup> Paradigm (multi-mode)
- SpectraMax<sup>®</sup> M5e/M5 (multi-mode)
- SpectraMax<sup>®</sup> M4/M3 (multi-mode)
- SpectraMax<sup>®</sup> M2e/M2 (dual-mode)
- Gemini<sup>™</sup> EM and Gemini<sup>™</sup> XPS (fluorescence)
- SpectraMax<sup>®</sup> L
- SpectraMax<sup>®</sup> Plus<sup>384</sup> (Serial#-MNR-xxxx)
- SpectraMax<sup>®</sup> 340PC<sup>384</sup> (Serial#-LNR-xxxx)
- SpectraMax<sup>®</sup> 190 (Serial#-NNR-xxxx)
- VersaMax<sup>™</sup> (Serial#-BNR-xxxx)

## Ordering information

Please contact your Molecular Devices representative for configuration options.

StakMax Microplate Stacker  
Part Number: STAKMAX

20-plate magazine (pair)  
Part Number: 0200-6016

40-plate magazine (pair)  
Part Number: 0200-6017

50-plate magazine  
Part Number: 0310-5102

Shortside barcode reader  
Part Number: 0310-5102

Longside barcode reader  
Part Number: 0310-5101

## Technical specifications

Performance specifications	
Compatibility	SBS-standard 96- and 384-well plates with plate height between 12–16 mm and 96-well "strip" plates
Barcode libraries	Interleaved 2 of 5, Code 128, EAN 128, Code 93, Code 39, Codabar
General specifications	
Dimensions (in.)	9.1 (H) x 17.3 (W) x 20.0 (D)
Dimensions (cm)	23.1 (H) x 44.0 (W) x 50.9 (D)
Bench space required (StakMax Stacker plus reader):	34.1 (W) x 20.7 (D) (in.)
	86.6 (W) x 52.6 (D) (cm)
Weight	26 lbs. (11.8 kg)
Power source	100–240 V <sub>AC</sub> , 3 A, 50/60 Hz, autoranging line voltage
Operating temp	10–40°C
Connectivity	USB 2.0 (requires SoftMax Pro Software v5.1 or greater for Windows)

## 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  
Brazil +55-11-3616-6607  
China (Beijing) +86-10-6410-8669  
China (Shanghai) +86-21-3372-1088  
Germany 00800-665-32860

Japan (Osaka) +81-6-7174-8831  
Japan (Tokyo) +81-3-6362-5260  
South Korea +82-2-3471-9531  
United Kingdom +44-118-944-8000

