

RapidStak automates SpectraMax microplate readers and AquaMax DW4 liquid handlers as stand-alone workstations

SPECTRAMAX TECH NOTE



While Molecular Devices and Thermo provide their respective products, the on-site integration and automation setup will be provided by the Thermo Scientific group.

PRODUCT INTEGRATION OPTIONS

From Molecular Devices

- SpectraMax absorbance readers
 - VersaMax™
 - SpectraMax 340PC³⁸⁴
 - SpectraMax 190
 - SpectraMax Plus³⁸⁴
- Gemini fluorometers
 - Gemini XPS
 - Gemini EM
- SpectraMax multi-mode readers
 - SpectraMax M2
 - SpectraMax M2^e
 - SpectraMax M5
 - SpectraMax M5^e
- SoftMax Pro 5 software
- AquaMax DW4 dispenser and washer

From Thermo Scientific

- RapidStak Microplate Stacker (30- or 50-plate capacity, up to 50 15-mm plate capacity)
- RapidStak Microplate Stacker 2x (up to 150 15-mm plate capacity)
- 25-, 30-, or 50-plate magazines available
- Polara RS hardware and software integration kit for Molecular Devices SpectraMax
- Polara RS hardware and software integration kit for Molecular Devices AquaMax DW4
- Polara RS hardware and software integration kit for Molecular Devices SpectraMax and AquaMax DW4

For more information please email info@moldev.com.

INTRODUCTION

The Thermo Scientific RapidStak® microplate stacker provides reliable walk-away automation to Molecular Devices' SpectraMax® microplate readers and AquaMax® DW4 liquid handlers. Users can run their ELISA or screening assays unattended with up to 150 15-mm capacity microplates.

Researchers can also integrate both a SpectraMax reader and an AquaMax DW4 liquid handler to a single RapidStak, enabling a dispense, wash, and read operation in a single workstation. The Polara RS automation software communicates with Molecular Devices SoftMax® Pro software enabling data acquisition, analysis and GxP capabilities and increasing the power of the workstation. The RapidStak features a unique choreography and telescoping arm design that allows it to deliver plates to the instrument while performing the next stack/de-stack operation simultaneously. This multi-tasking yields a load/unload time of less than six seconds.

SALES AND SUPPORT

United States & Canada
Molecular Devices
Tel. +1-800-635-5577
Fax +1-408-747-3601

Australia
Molecular Devices Pty. Ltd.
Tel. +61-3-9896-4700
Fax +61-3-8640-0742

Brazil
Molecular Devices Brazil
Tel. +55-11-3616-6607
Fax +55-11-3616-6607

China
Molecular Devices Beijing
Tel. +86-10-6410-8669
Fax +86-10-6410-8601

Molecular Devices Hong Kong
Tel. +85-2-8125-2509
Fax +85-2-3010-2828

Molecular Devices Shanghai
Tel. +86-21-6887-8820
Fax +86-21-6887-8890

Germany
Molecular Devices GmbH
Tel. +49-89-9605-880
Fax +49-89-9620-2345

Japan
Nihon Molecular Devices Osaka
Tel. +81-6-6399-8211
Fax +81-6-6399-8212

Nihon Molecular Devices Tokyo
Tel. +81-3-5282-5261
Fax +81-3-5282-5262

South Korea
Molecular Devices Korea, LLC
Tel. +82-2-3471-9531
Fax +82-2-3471-9532

United Kingdom
Molecular Devices Ltd.
Tel. +44-118-944-8000
Fax +44-118-944-8001

www.moleculardevices.com

AquaMax, Molecular Devices, SoftMax, and SpectraMax are registered trademarks of MDS Analytical Technologies. All other trademarks are the property of their respective owners.
©2007 MDS Analytical Technologies. Printed in U.S.A.
5/07 EP #0120-1370A

