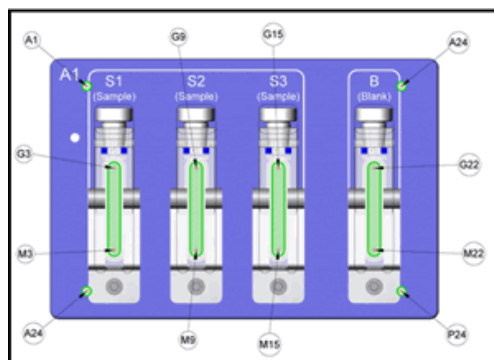


SpectraCuvette Adapter

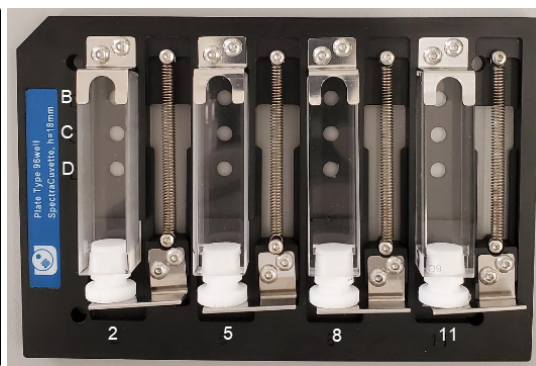
The SpectraCuvette Adapters from Molecular Devices are designed to allow microplate readers to read cuvettes.

Cuvette Part Numbers and Tested Instruments

Part Number	Height	Well Design	Tested Microplate Readers
0200-7204	22 mm	384	FilterMax™ F3 & F5 Multi-Mode Microplate Readers, SpectraMax® i3x Multi-Mode Detection Platform, SpectraMax® iD3 & iD5 Multi-Mode Microplate Readers, and SpectraMax® Paradigm® Multi-Mode Microplate Reader
5301193	18 mm	96	SpectraMax i3x, SpectraMax iD3, SpectraMax iD5, SpectraMax M Series, and SpectraMax® Mini Multi-Mode Microplate Reader



0200-7204 384-well



5301193 96-well

Read Areas for the Cuvettes

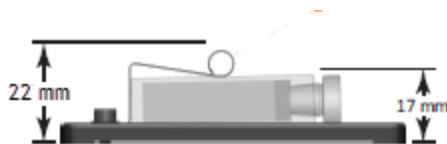
Each cuvette adapter has positions marked for samples and/or a blank.

Part Number 0200-7204		Part Number 5301193	
Cuvette Position	Read Area 384-well plate	Cuvette Position	Read Area 96-well plate
Sample 1 (S1)	Column 3 (G3 through M3)	Sample 1 (S1)	Column 2 (B2 through D2)
Sample 2 (S2)	Column 9 (G9 through M9)	Sample 2 (S2)	Column 5 (B5 through D5)
Sample 3 (S3)	Column 15 (G15 through M15)	Sample 3 (S3)	Column 8 (B8 through D8)
Blank (B)	Column 22 (G22 through M22)	Sample 4 (S4)	Column 11 (B11 through D11)

The notched corner indicates the A1 corner, and alignment holes work with the Microplate Optimization Wizard in the SoftMax® Pro Data Acquisition and Analysis Software to accurately align the instrument with the adapter. Place the adapter onto the plate drawer with the label side up.

Read Height Clearance

Older 384-well SpectraCuvette Adapters (part number 0200-7204 without a silver sticker) have a plate height of 24 mm and you cannot use this in the SpectraMax iD3 or SpectraMax iD5 instrument because these instruments have a maximum plate height of 22 mm. To prevent damage to the instrument, use the correct adapter (22 mm or 18 mm) as listed in the table on the previous page.



Insert Cuvettes into the SpectraCuvette Adapter

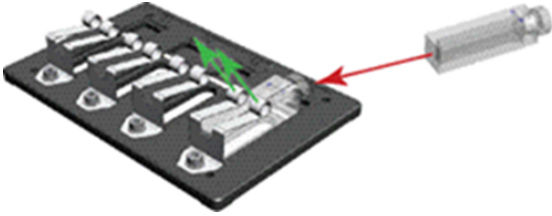
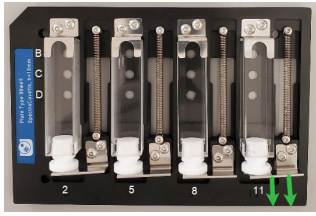
You can insert up to four cuvettes in the SpectraCuvette Adapter. The cuvettes must be 12.5 mm W x 12.5 mm D x ~46 mm H and have a stopper.



CAUTION! Make sure to properly close the cuvette stopper to prevent leakage. Do not let water or other fluids drip inside the instrument.



WARNING! PINCH HAZARD. Do not place your finger between the retaining spring and the cuvette.

384-well plate 0200-7204	96-well plate 5301193
	
<p>1. Gently pull up the retaining spring where you want to insert the cuvette.</p>	<p>1. Gently pull down on the retaining spring where you want to insert the cuvette.</p>
<p>2. Place the cuvette completely into the recessed area with the BOTTOM of the cuvette against the bottom of the retaining spring.</p>	<p>2. Place the cuvette completely into the recessed area with the TOP of the cuvette against the retaining spring.</p>
<p>3. Gently lower the retaining spring to secure the cuvette into position.</p>	<p>3. Gently release retaining spring to secure the cuvette into position.</p>

For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/patents>

Product manufactured by Molecular Devices, LLC.

3860 N. First Street, San Jose, California, 95134, United States of America.

Molecular Devices, LLC is ISO 9001 registered.

©2022 Molecular Devices, LLC.

All rights reserved.