

Unlocking the SpectraMax Paradigm Instrument

When the SpectraMax® Paradigm® Multi-Mode Detection Platform is first installed, the hardware transport locks need to be removed. As an additional safety precaution, internal locks controlled by the software prevent the drawers from opening until the instrument is detected and initialized by the controlling software.



Note: Do not install detection cartridges until after you have completed the following instructions to unlock the drawers and initialize the instrument.

To unlock the drawers and initialize the instrument:

1. Make sure that the hardware transport locks have been removed. See the SpectraMax Paradigm Instrument unpacking guide or user guide.

CAUTION! The instrument can be damaged if all the transport locks are not removed before unlocking the drawers and initializing the instrument.

2. Make sure that all drawers have been pushed back into the instrument, the front panel has been replaced, and the microplate door is closed.
3. Make sure that the instrument is connected to the host computer and to a power source, and that the SoftMax Pro Software has been installed on the host computer. See the SpectraMax Paradigm Instrument unpacking guide or user guide.
4. Turn on the power switch on the back of the instrument.

If the **Standby** button on the front lower-right corner of the instrument is illuminated, press the button to take the instrument out of standby mode.

The LEDs on the status panel flash and turn off, and then the amber LED turns on indicating that the instrument drawers are locked.

5. Start the SoftMax Pro Software as described in the *SoftMax Pro Software User Guide*.

The amber LED on the status panel turns off and the green LED turns on, indicating a successful connection between the instrument and the software.

6. After successfully connecting to the instrument, the **Instrument Unlocking Procedure** wizard appears.

If the **Instrument Unlocking Procedure** wizard does not appear, then use the **Instrument Connection** dialog to select and connect to the instrument as described in the *SoftMax Pro Software User Guide*.



Note: If the instrument does not appear in the **Available Instruments** list of the **Instrument Connection** dialog, then click **Refresh** above the list.

7. Follow the on-screen instructions in the **Instrument Unlocking Procedure** wizard to unlock the drawers and initialize the instrument.

After completing the **Instrument Unlocking Procedure** wizard, the instrument performs an initialization procedure that moves the optics and microplate drawers to home positions. The green LED on the status panel turns on, the amber LED flashes and turns off, and then the green LED remains on.

This document is provided to customers who have purchased Molecular Devices, LLC ("Molecular Devices") equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a license agreement. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Molecular Devices makes no warranties or representations as to the fitness of this equipment for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom.



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/productpatents>

Product manufactured by Molecular Devices, LLC.
1311 Orleans Drive, Sunnyvale, California, United States of America 94089.
Molecular Devices, LLC is ISO 9001 registered.
© 2012 Molecular Devices, LLC.
All rights reserved.
