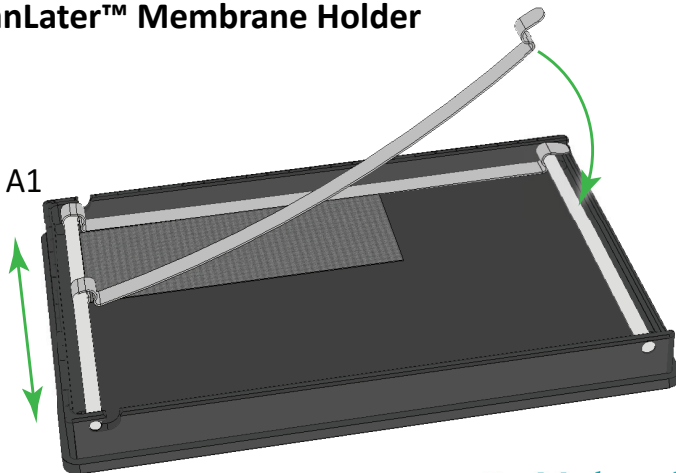


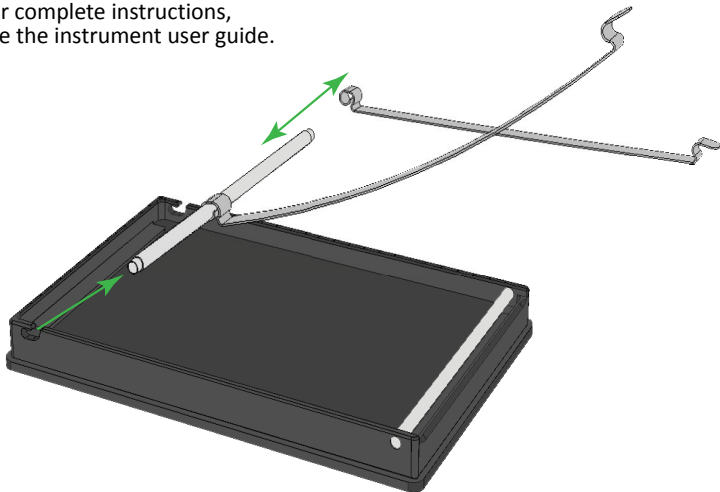
ScanLater™ Membrane Holder



ED 000 104

 **Molecular
Devices**

For complete instructions,
see the instrument user guide.



© 2013 Molecular Devices, LLC. All rights reserved.
For research use only. Not for use in diagnostic procedures.