



## Certificate of Compliance

Issue Date: February 5, 2020

Molecular Devices, LLC certifies that SoftMax<sup>®</sup> Pro, SoftMax<sup>®</sup> Pro GxP, and GxP Admin software (“Software”) were tested and verified to be in compliance with all of our internal, applicable manufacturing and quality requirements.

Molecular Devices, LLC certifies that the Software was designed, developed, and verified per formal documented processes to meet all functional and performance specifications. To support this declaration, all the supporting documents are archived and maintained by Molecular Devices, LLC and are available for reference and review at our corporate offices in San Jose, California, U.S.A.

SoftMax<sup>®</sup> Pro GxP software was designed to be used as part of a 21 CFR Part 11 compliant system. The compliance relies on an administrator’s ability to create a secure environment for generating, analyzing, and storing data.

The above systems, processes and documentation are formal requirements of the Molecular Devices, LLC Quality Management System.



ISO 9001  
FS 534246

Signed:

A handwritten signature in blue ink, appearing to read "Shreen Ibrahim", written over a horizontal line.

Shreen Ibrahim  
Director, Global Quality and EHS  
Molecular Devices, LLC  
3860 N First Street  
San Jose, CA 95134