



## DECLARATION OF CONFORMITY

(According to Decision 768/2008)



**Manufacturer's Name:** Molecular Devices, LLC

**Manufacturer's Address:** 3860 N. 1<sup>st</sup> Street, San Jose, CA 95134 USA

**This declaration of conformity is issued under the sole responsibility of the manufacturer.**

**Product Name:** Spectrophotometer

**Model Numbers:** SpectraMax QuickDrop

**Product Options:** This declaration covers all options of the above product(s)

**The object of the declaration described above is in conformity with:**

Low Voltage Directive (2014/35/EU)  
EMC Directive (2014/30/EU)  
RoHS Directive (2011/65/EU)  
REACH Regulation (EC No 1907/2006)

**Harmonized standards used:**

**EMC** IEC 61326-1:2012 / EN 61326-1:2013

**LVD** IEC 61010-1:2010 / EN 61010-1:2010

**RoHS** EN 50581:2012

I, the undersigned, hereby declare that the equipment specified above conforms to the above Directives and Standards.

March 11, 2021

Date

San Jose, CA

Place of Signature

Director, Global Quality