



DECLARATION OF CONFORMITY

(According to Decision 768/2008)



Manufacturer's Name: Molecular Devices, LLC

Manufacturer's Address: 3860 N. First Street, San Jose, CA 95134, USA

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

Product Name: Single Cell Printer

Model Numbers: Cloneselect Single-Cell Printer

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)

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.

July 18, 2018

Date

Sunnyvale, CA

Place of Signature

Director, Global Quality