

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



EU Representative: Molecular Devices (Austria) GmbH
EU Representative Address: 17 Urstein Süd, Puch / Hallein Salzburg, Austria 5412

UK Representative: Molecular Devices (UK) Limited
UK Representative Address: 660-665 Eskdale Road, Winnersh Triangle, Wokingham, Berkshire RG41 5TS, United Kingdom

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

Product Name: MicroArray System
Model Numbers: GenePix Series
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 (EU) 2015/863
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 IEC 63000:2018/EN 63000:2018

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

May 9, 2022

Date

San Jose, CA

Place of Signature



Director, Global Quality