

DECLARATION OF CONFORMITY

(According to Decision 768/2008)

| | | |
|--|---|---|
| <u>Manufacturer's Name:</u> | Molecular Devices, LLC |  |
| <u>Manufacturer's Address:</u> | 3860 N. First Street, San Jose, CA 95134, USA | |
| <u>EU Representative:</u> | Molecular Devices (Austria) GmbH | |
| <u>EU Representative Address:</u> | 17 Urstein Süd, Puch / Hallein Salzburg, Austria 5412 | |
| <u>UK Representative:</u> | Molecular Devices (UK) Limited | |
| <u>UK Representative Address:</u> | 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.

| | |
|--------------------------------|---|
| <u>Product Name:</u> | Microplate Reader |
| <u>Model Numbers:</u> | SpectraMax iD3, iD5 Series |
| <u>Product Options:</u> | 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)
Machine Directive (2006/42/EC)
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 61010-1:2010/61010-2-081 EN 61010-1:2010/EN 61010-2-081 |
| 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