

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:

1180 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:

Multi-Mode Microplate Reader

Model Numbers:

SpectraMax i3x w/MiniMax 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)

Machine Directive (2006/42/EC)

EMC Directive (2014/30/EU)

Packaging and Packaging Waste Regulation 2025/40 (PPWR)

RoHS Directive (EU) 2015/863

REACH Regulation (EC No 1907/2006)

Harmonized standards used:

EMC IEC/EN 61326-1:2021

LVD IEC/EN 61010-1/61010-2-081+A1/60825-1

RoHS IEC/EN 63000:2018

Packaging CR 13695-1:2020

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

July 1, 2026

Date

San Jose, CA

Place of Signature

Director, Global Quality