


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>	1180 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>	Hypoxic Chamber Bacteria Picker
<u>Model Numbers:</u>	QPix FLEX 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/EN 61326-1:2021
LVD	IEC 61010-1:2010/AMD1:2016
RoHS	IEC/EN 63000:2018

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

April 21, 2026
Date

San Jose, CA
Place of Signature



Director, Global Quality