


## **DECLARATION OF CONFORMITY**

(According to Decision 768/2008)

<b><u>Manufacturer's Name:</u></b>	Molecular Devices, LLC	
<b><u>Manufacturer's Address:</u></b>	3860 N. First Street, San Jose, CA 95134, USA	
<b><u>EU Representative:</u></b>	Molecular Devices (Austria) GmbH	
<b><u>EU Representative Address:</u></b>	17 Urstein Süd, Puch / Hallein Salzburg, Austria 5412	
<b><u>UK Representative:</u></b>	Molecular Devices (UK) Limited	
<b><u>UK Representative Address:</u></b>	1180 Eskdale Road, Winnersh Triangle, Wokingham, Berkshire RG41 5TS, United Kingdom	

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

<b><u>Product Name:</u></b>	Automated Cell Culture System
<b><u>Model Numbers:</u></b>	CellXpress.ai
<b><u>Product Options:</u></b>	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:**

<b>EMC</b>	IEC 61326-1:2020/EN 61326-1:2021
<b>LVD</b>	IEC 61010-1:2010/AMD1:2016
<b>RoHS</b>	IEC 63000:2018/EN 63000:2018

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

February 27, 2026

Date

San Jose, CA

Place of Signature



Director, Global Quality