

EC Declaration of Conformity

Manufacturer Name:	Cyvek, Inc.	
Manufacturer Address:	5 Technology Drive Wallingford, CT 06492 USA	
Manufacturer SRN: (Single Registration Number)	US-MF-000049483	
EU Authorized Representative:	MedEnvoy Global BV Prinses Margrietplantsoen 33, Suite 123 2595 AM The Hague The Netherlands	
Product Name:	Ella	
Product Catalogue Number(s) / Code(s)	<table border="1"><tr><td>REF</td></tr></table> 600-100-CE	REF
REF		
Basic UDI-DI	86001420031000139001CEYD	
Classification	Class A non-sterile, Rule 5(b)	
Intended Purpose	<p>The Ella Instrument is a non-automated instrument intended for in vitro diagnostic use with a supported Ella IVD Assay to perform immunoassay diagnostics.</p> <p>The Ella Instrument sequences the processing of samples, reagents and wash buffers, and measures and sorts fluorescence signals, using a supported Ella IVD Assay Kit.</p> <p>The Ella Instrument is intended for use by trained laboratory personnel.</p>	

Common Specifications / Standards Applied ISO 13485:2016
IEC 61010-1:2010 / AMD1:2016
IEC 61326:2021
EN ISO 14971:2019
IEC 60825-1:2014
IEC 62304:2006/A1:2015
IEC 62366-1:2015

Directives EC 1907 / 2006
2011 / 65 / EU
2015 / 863 / EU

This declaration of conformity is issued under the sole responsibility of Cyvek, Inc. We hereby declare that the medical device(s) specified above meet the provision of the Regulation (EU) IVDR 2017/746 for medical devices. This declaration is supported by the Cyvek Inc. Quality Management System's certification to ISO 13485:2016.

All supporting documentation is retained by the manufacturer.

Roberto L. Muñoz

Roberto Muñoz
Director, Quality Assurance

Wallingford, Connecticut, U.S.A
January 22, 2026

Place and date of Issue