



Meso Scale Diagnostics, LLC.

PACKAGING DECLARATION OF CONFORMITY

MSD® # D0000-6523-A

This Declaration of Conformity is issued under the sole responsibility of the manufacturer. The packaging products listed are in conformity with the applicable Union harmonization legislation: Packaging and Packaging Waste Regulation (PPWR) (EU) 2025/40 as amended and its applicable requirements.

Manufacturer Name: Meso Scale Diagnostics, LLC.
Manufacturer's Address: 1601 Research Boulevard
Rockville, Maryland 20850 USA

Declares that the packaging families and covered items:

Packaging Family Name: Cardboard and corrugate
Packaging Description: Solid cardboard, corrugated cardboard, kraft cardboard instrument and consumable packaging

Conforms to the following Union harmonization legislation:

The requirements specified in Annex VII of PPWR for maintaining technical documentation have been met.

The requirements specified in Article 5 of PPWR for verification that the cumulative concentration levels of Lead, Cadmium, Mercury, and Hexavalent Chromium present in the packaging do not exceed 100 mg/kg.

The requirements specified in Article 44 and 45 of PPWR requiring producers to register with the national registers and report packaging volume data for each packaging type have been met.

As manufacturer we declare that the product packaging follows the provisions of the standards and specifications stated above. We are actively monitoring upcoming PPWR deadlines to ensure compliance as implementation dates are announced.

EAR Name: MedEnvoy Global B.V
EAR Address: Prinses Margrietplantsoen 33, Suite 123
2595 AM The Hague, The Netherlands

Date of Issue: 04 AUG 2026
Place of Issue: Rockville, MD, USA

Lillian J. Quintero
VP, Compliance & Bioanalytical Testing