

**CERTIFICATE OF CONFORMANCE**

Product Description: Mouse Isotyping Panel 1 Kit
Catalog Number: K15183B-Series
Lot Number: K0050937
Expiration Date See expiration date on specific components below

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse Isotyping Panel 1 96-Well 10-Spot SECTOR Plate	Z0057677	2-8°C	30 SEP 2028
Mouse IgA Antibody	D0050916	2-8°C	30 SEP 2029
Mouse IgG1 Antibody	D0050917	2-8°C	30 SEP 2029
Mouse IgG2a Antibody	D0050918	2-8°C	30 SEP 2029
Mouse IgG2b Antibody	D0050919	2-8°C	31 OCT 2029
Mouse IgG3 Antibody	D0050920	2-8°C	30 SEP 2029
Mouse IgM Antibody	D0050921	2-8°C	30 SEP 2029
Mouse Isotyping Panel 1 Calibrator Blend	A0020309	≤ -70°C	31 MAY 2028

Plate Specifications:

Coating Uniformity	Specifications
Mean Intraplate CV	≤ 15%
Mean Interplate CV	≤ 25%
No plate with CV	> 18%

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 JUN 2026