



CERTIFICATE OF ANALYSIS

COA - 07179

Product Description: MSD[®] Mouse Isotyping Panel 1 Kit

Catalog #/Size: K15183B-1 / 1 Plate; K15183B-2 / 5 Plates; K15183B-4 / 25 Plates

Lot Number: K0050662

Expiration Date: N/A (See expiration date on specific components below)

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Kit Contents:

Plate Lot Number: Z0054938

Plate Lot Catalog Number: N75183A

Plate Lot Expiration Date: 31 AUG 2016

Antibody blend Lot Number: D0050570

Antibody blend Catalog Number: D22JJ

Antibody blend Expiration Date: 31 JUL 2016

Antibody blend Lot Number: D0050571

Antibody blend Catalog Number: D22JL

Antibody blend Expiration Date: 31 JUL 2016

Antibody blend Lot Number: D0050572

Antibody blend Catalog Number: D22JM

Antibody blend Expiration Date: 31 JUL 2016

Antibody blend Lot Number: D0050573

Antibody blend Catalog Number: D22JN

Antibody blend Expiration Date: 31 AUG 2016

Antibody blend Lot Number: D0050574

Antibody blend Catalog Number: D22JO

Antibody blend Expiration Date: 31 JUL 2016

Antibody blend Lot Number: D0050575

Antibody blend Catalog Number: D22JP

Antibody blend Expiration Date: 31 JUL 2016

Calibrator blend Lot Number: A0020156

Calibrator blend Catalog Number: C0183

Calibrator blend Expiration Date: 30 SEP 2016



CERTIFICATE OF ANALYSIS

COA - 07179

Specifications:

Release method: Coating uniformity

Mean Intraplate CV: $\leq 15\%$

Mean Interplate CV: $\leq 18\%$

No more than one plate with CV $> 18\%$

No plate with CV $> 20\%$

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Xiomara Talavera	Quality Control	<i>Talavera</i>	02 AUG 2016
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	03 AUG 2016