



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD® Mouse Isotyping Panel 1 Kit

**Catalog Number:** K15183B - Series

**Lot Number:** K0050835

**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					
<b>Lot #</b>	Z0056321	<b>Catalog #</b>	N75183A-1	<b>Expiration Date:</b>	28 FEB 2021

Antibody Blend	
<b>Lot Number:</b>	D0050726
<b>Catalog Number:</b>	D22JJ-2
<b>Expiration Date:</b>	29 FEB 2020
<b>Storage Temperature:</b>	2-8°C

Antibody Blend	
<b>Lot Number:</b>	D0050730
<b>Catalog Number:</b>	D22JO-2
<b>Expiration Date:</b>	29 FEB 2020
<b>Storage Temperature:</b>	2-8°C

Antibody Blend	
<b>Lot Number:</b>	D0050727
<b>Catalog Number:</b>	D22JL-2
<b>Expiration Date:</b>	30 APR 2020
<b>Storage Temperature:</b>	2-8°C

Antibody Blend	
<b>Lot Number:</b>	D0050731
<b>Catalog Number:</b>	D22JP-2
<b>Expiration Date:</b>	29 FEB 2020
<b>Storage Temperature:</b>	2-8°C

Antibody Blend	
<b>Lot Number:</b>	D0050728
<b>Catalog Number:</b>	D22JM-2
<b>Expiration Date:</b>	29 FEB 2020
<b>Storage Temperature:</b>	2-8°C

Calibrator Blend	
<b>Lot Number:</b>	A0020225
<b>Catalog Number:</b>	C0183-2
<b>Expiration Date:</b>	29 FEB 2020
<b>Storage Temperature:</b>	≤ -70°C

Antibody Blend	
<b>Lot Number:</b>	D0050729
<b>Catalog Number:</b>	D22JN-2
<b>Expiration Date:</b>	29 FEB 2020
<b>Storage Temperature:</b>	2-8°C

Specifications:

Release method:

Coating Uniformity	Specifications
Mean Intraplate CV	≤ 15%
Mean Interplate CV	≤ 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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	16 NOV 2018