



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD<sup>®</sup> Mouse MIP-3 $\alpha$  Kit

**Catalog Number:** K152MSD-Series

**Lot Number:** K00E0261

**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 #	Z00E0211	Catalog #	N452MSA	Expiration Date:	31 JUL 2018

Antibody Blend	
LotNumber:	D00E0241
Catalog Number:	D22MS
Expiration Date:	31 AUG 2018
Storage Temperature:	2-8°C

Calibrator Blend	
Lot Number:	A00E130
Catalog Number:	C02MS
Expiration Date:	31 AUG 2018
Storage Temperature:	$\leq -70^{\circ}\text{C}$

Specifications:

Release method:

Coating Uniformity	Specifications
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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	13 APR 2017