



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Mouse GM-CSF Ultra-Sensitive Kit

Catalog Number: K152AXC-1 / 1 Plate; K152AXC-2 / 5 Plates; K152AXC-4 / 25 Plates

Lot Number: K0035187

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	Z0046128	Catalog #	N45022B	Expiration Date	31 OCT 2018

Antibody Blend					
Lot	D0033652	Catalog #	D22AE	Expiration Date	31 OCT 2018
Spot #	Description		Host / Species		
1	mIFNg		Rat		

Calibrator Blend					
Lot	A0032245	Catalog #	C0022	Expiration Date	30 NOV 2018
Spot #	Description		Host / Species		
1	Mouse GM-CSF		<i>E. coli</i>		

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
Preparer	Xiomara Talavera	Quality Control	<i>X Talavera</i>	05 DEC 2016
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	06 DEC 2016