



CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: MSD® Mouse ProInflammatory 7-Plex Tissue Culture Kit

Catalog Number: K15012B - Series

Lot Number: K0035288

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	Z0046254	Catalog #	N75012B	Expiration Date	31 MAR 2019

Antibody Blend					
Lot	D0033728	Catalog #	D2012	Expiration Date	31 MAR 2019
Spot #	Description		Host / Species		
1	mIL-1beta		Goat		
2	mIL-12		Goat		
3	Mouse IFN-gamma		Rat		
4	Mouse IL-6		Goat		
5	Mouse KC/GRO		Goat		
6	Mouse IL-10		Goat		
7	mTNFa		Rat		

Calibrator Blend					
Lot	A0032278	Catalog #	C0012	Expiration Date	31 MAR 2019
Spot #	Description		Host / Species		
1	Mouse IFN-g		<i>E. coli</i>		
2	Mouse IL-1b		<i>E. coli</i>		
3	Mouse IL-6		<i>E. coli</i>		
4	Mouse KC/GRO		<i>E. coli</i>		
5	Mouse IL-10		<i>E. coli</i>		
6	Mouse IL-12t/p70		<i>Sf 21</i>		
7	Mouse TNF-a		<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%

COA-07856-A



CERTIFICATE OF ANALYSIS

Page 2 of 2

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>	06 FEB 2017
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	07 FEB 2017