



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K15012B - Series

Lot Number: K0035338

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 #	Z0046333	Catalog #	N75012B	Expiration Date:	31 JUL 2019

Antibody Blend	
Lot Number:	D0033761
Catalog Number:	D2012
Expiration Date:	30 JUN 2019
Storage Temperature:	2-8°C
Spot #	Description
1	mIL-1beta
2	mIL-12
3	Mouse IFN-gamma
4	Mouse IL-6
5	Mouse KC/GRO
6	Mouse IL-10
7	mTNFa

Calibrator Blend	
Lot Number:	A0032254
Catalog Number:	C0012
Expiration Date:	31 DEC 2018
Storage Temperature:	≤ -70°C
Spot #	Description
1	Mouse ProInflammatory Panel

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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	25 JAN 2018