



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human ProInflammatory I 4 Plex Tissue Culture Kit

Catalog Number: K15009B - Series

Lot Number: K0035333

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 #	Z0046320	Catalog #	N45009B	Expiration Date:	31 MAY 2019

Antibody Blend	
Lot Number:	D0033758
Catalog Number:	D2009
Expiration Date:	31 MAY 2019
Storage Temperature:	2-8°C
Spot #	Description
1	Human IFN-gamma
2	Human IL-1β
3	Human IL-6
4	Human TNF-alpha

Calibrator Blend	
Lot Number:	A0032258
Catalog Number:	C0105
Expiration Date:	31 JAN 2019
Storage Temperature:	≤ -70°C
Spot #	Description
1	Human Tissue Culture Calibrator 2

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>	08 NOV 2017