



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Mouse TH1/TH2 9-Plex Tissue Culture Kit

Catalog Number: K15013B - Series

Lot Number: K0035444

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 #	Z0046457	Catalog #	N05013B	Expiration Date:	30 SEP 2019

Antibody Blend	
Lot Number:	D0033842
Catalog Number:	D2013
Expiration Date:	30 NOV 2019
Storage Temperature:	2-8°C
Spot #	Description
1	mIFN-gamma
2	mIL-1beta
3	mIL-2
4	mIL-4
5	m/rIL-5
6	mKC/GRO/CINC/CXCL1
7	mIL-10
8	mIL-12
9	mTNF-alpha

Calibrator Blend	
Lot Number:	A0032314
Catalog Number:	C0013
Expiration Date:	30 SEP 2019
Storage Temperature:	≤ -70°C
Spot #	Description
1	Mouse TH1/TH2 9-Plex

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>	18 JUN 2018