



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K15013B - Series

Lot Number: K0034851

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 #	Z0045597	Catalog #	N05013B	Expiration Date:	31 JUL 2017

Antibody Blend	
Lot Number:	D0033365
Catalog Number:	D2013
Expiration Date:	31 JUL 2017
Storage Temperature:	2-8°C
Spot #	Description
1	Mouse IFN-gamma
2	mIL-1beta
3	Mouse IL-2
4	Mouse IL-4
5	m/r IL-5
6	Mouse KC/GRO
7	mIL-10
8	Mouse IL-12
9	mTNFa

Calibrator Blend	
Lot Number:	A0032111
Catalog Number:	C0013
Expiration Date:	31 JUL 2017
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	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	25APR2017