



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human TH1/TH2 10-Plex Tissue Culture Kit

Catalog Number: K15010B-Series

Lot Number: K0035328

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 #	Z0046347	Catalog #	N05010B	Expiration Date:	30 JUN 2019

Antibody Blend	
Lot Number:	D0033765
Catalog Number:	D2010
Expiration Date:	30 JUN 2019
Storage Temperature:	2-8°C
Spot #	Description
1	Human TNF-alpha
2	hIFN-gamma
3	Human IL-1beta
4	Human IL-2
5	hIL-4
6	Human IL-5
7	Human IL-8
8	hIL-10
9	hIL12p70
10	hIL-13

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	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	10MAY2017