



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human TNF- α Tissue Culture Kit

Catalog Number: K151BHB-Series

Lot Number: K0035293

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 #	Z0046266	Catalog #	L451BHB	Expiration Date:	30 APR 2019

Antibody Blend	
Lot Number:	D0033733
Catalog Number:	D21BH
Expiration Date:	31 MAR 2019
Storage Temperature:	2-8°C
Spot #	Description
1	Human TNF α

Calibrator Blend	
Lot Number:	A0032258
Catalog Number:	C0105
Expiration Date:	31 JAN 2019
Storage Temperature:	$\leq -70^{\circ}\text{C}$
Spot #	Description
1	Human Tissue Culture Calibrator 2

Specifications:

Release method:

Coating Uniformity	Specifications
Mean Intraplate CV	$\leq 15\%$
Mean Interplate CV	$\leq 18\%$
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>	13 APR 2017