



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human TNF-RI Ultra-Sensitive Kit

Catalog Number: K151BIC-Series

Lot Number: K0035531

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 #	Z0046594	Catalog #	N45016A	Expiration Date:	31 JAN 2020

Antibody Blend	
Lot Number:	D0033861
Catalog Number:	D21B1
Expiration Date:	30 NOV 2019
Storage Temperature:	2-8°C
Spot #	Description
1	Human TNF-RI

Calibrator Blend	
Lot Number:	A0032341
Catalog Number:	C0016
Expiration Date:	31 DEC 2019
Storage Temperature:	≤ -70°C
Spot #	Description
1	hTNF-RII

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>	13 DEC 2017