



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human IFN-γ Tissue Culture Kit

Catalog Number: K151AEB-Series

Lot Number: K0035619

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 #	Z0046678	Catalog #	L451AEB	Expiration Date:	31 MAR 2020

Antibody Blend	
LotNumber:	D0033974
Catalog Number:	D21AE
Expiration Date:	31 MAR 2020
Storage Temperature:	2-8°C
Spot #	Description
1	Human IFN γ

Calibrator Blend	
Lot Number:	A0032308
Catalog Number:	C01AE
Expiration Date:	30 SEP 2019
Storage Temperature:	$\leq -70^{\circ}\text{C}$
Spot #	Description
1	hIFN-g

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>	27 AUG 2018