



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® Human IFN-alpha2a Ultra-Sensitive Kit

Catalog Number: K151ACC-Series

Lot Number: K0034661

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	Z0045351	Catalog #	N45032A	Expiration Date	31 DEC 2016

Antibody Blend					
Lot	D0033231	Catalog #	D21AC	Expiration Date	31 DEC 2016
Spot #	Description		Host / Species		
1	h IFN-alpha2a		Mouse		

Calibrator Blend					
Lot	A0032044	Catalog #	C0032	Expiration Date	31 DEC 2016
Spot #	Description		Host / Species		
1	hIFN-a		<i>E. coli</i>		

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
Preparer	Xiomara Talavera	Quality Control	<i>Talavera</i>	08 DEC 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	12 DEC 2016