



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** K151ACC-Series

**Lot Number:** K0034869

**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					
<b>Lot</b>	Z0045626	<b>Catalog #</b>	N45032A	<b>Expiration Date</b>	31 AUG 2017

Antibody Blend					
<b>Lot</b>	D0033383	<b>Catalog #</b>	D21AC	<b>Expiration Date</b>	31 JUL 2017
<b>Spot #</b>	<b>Description</b>		<b>Host / Species</b>		
1	h IFN-alpha2a		Mouse		

Calibrator Blend					
<b>Lot</b>	A0032118	<b>Catalog #</b>	C0032	<b>Expiration Date</b>	31 JUL 2017
<b>Spot #</b>	<b>Description</b>		<b>Host / Species</b>		
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
<b>Preparer</b>	Xiomara Talavera	Quality Control	<i>X Talavera</i>	16 JAN 2017
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	17 JAN 2017