

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** U-PLEX® and U-PLEX Plus Human  $\alpha$ -Synuclein Kit  
**Kit Catalog Numbers:** K151WK Series  
**Kit Lot Number:** K0080739  
**Expiration Date:** 28 Feb 2019

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Human $\alpha$ -Synuclein Control 1		Human $\alpha$ -Synuclein Control 2		Human $\alpha$ -Synuclein Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Human $\alpha$ -Synuclein	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Human $\alpha$ -Synuclein	8.00	80% - 120%	< 20%	6800	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Human $\alpha$ -Synuclein	20X Calibrator	210000	pg/mL
		Human $\alpha$ -Synuclein Control 1	17500	pg/mL
		Human $\alpha$ -Synuclein Control 2	1710	pg/mL
		Human $\alpha$ -Synuclein Control 3	209	pg/mL

Controls tested may not be the same lot as shipped.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

### Standard Curve Data:

Human $\alpha$ -Synuclein - Spot 1		
Conc. pg/mL	ECL Signal	%CV
10500	456,000	2.9
2630	96,300	3.6
656	20,900	3.4
164	4,920	6.2
41.0	1,270	5.8
10.3	371	5.5
2.56	143	5.4
0.00	67	5.9

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD Gold 96-Well Small Spot Streptavidin Plate	Z0021268	2-8°C	31 Jul 2019
Biotin Anti-hu $\alpha$ -Synuclein	D0080944	2-8°C	28 Feb 2019
SULFO-TAG Anti-hu $\alpha$ -Synuclein	D0080945	2-8°C	31 Aug 2019
Human $\alpha$ -Synuclein Calibrator	A0080115	$\leq -70^{\circ}\text{C}$	30 Jun 2019
Diluent 49	M0490003	$\leq -10^{\circ}\text{C}$	30 Apr 2019
Diluent 49	M0490003	$\leq -10^{\circ}\text{C}$	30 Apr 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in U-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Human $\alpha$ -Synuclein Control 1	Not Kit Specific
Human $\alpha$ -Synuclein Control 2	Not Kit Specific
Human $\alpha$ -Synuclein Control 3	Not Kit Specific

### Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	27 FEB 2019