

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: U-PLEX® and U-PLEX Plus Human α -Synuclein Kit
Kit Catalog Numbers: K151WK Series
Kit Lot Number: K0080765
Expiration Date: 30 Jun 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Human α -Synuclein Control 1		Human α -Synuclein Control 2		Human α -Synuclein Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Hu α -Synuclein	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Hu α -Synuclein	8.00	80% - 120%	< 20%	6800	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Hu α -Synuclein	20X Calibrator	198200	pg/mL
		Human α -Synuclein Control 1	17923	pg/mL
		Human α -Synuclein Control 2	1690	pg/mL
		Human α -Synuclein Control 3	178	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:

Hu α -Synuclein - Spot 1		
Conc. pg/mL	ECL Signal	%CV
9910	539,000	3.1
2480	118,000	2.7
619	24,700	2.6
155	5,580	11.0
38.7	1,490	2.8
9.68	447	4.4
2.42	182	5.0
0.00	96	5.4

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD Gold 96-Well Small Spot Streptavidin Plate	Z0021280	2-8°C	31 Aug 2019
Biotin Anti-Hu α -Synuclein Ab	D0080978	2-8°C	30 Jun 2019
SULFO-TAG Anti-Hu α -Synuclein Ab	D0080979	2-8°C	30 Nov 2019
Human α -Synuclein Calibrator	A0080128	\leq -70°C	30 Sep 2019
Diluent 49	M0490004	\leq -10°C	30 Sep 2019
Diluent 49	M0490004	\leq -10°C	30 Sep 2019
MSD Read Buffer T (4X)	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 α -Synuclein Control 1	Not Kit Specific
Human α -Synuclein Control 2	Not Kit Specific
Human α -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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	27 FEB 2019