

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: K0081002
Expiration Date: 30 Sep 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	208000	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
10400	461,000	6.3
2600	94,400	4.2
650	19,700	3.7
163	4,720	4.5
40.6	1,250	3.2
10.2	372	4.1
2.54	157	2.9
0.00	90	9.8

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD Gold 96-well Small Spot Streptavidin SECTOR Plate	Z0021361	2-8°C	30 Apr 2020
SULFO-TAG Anti-hu α -Synuclein	D0081085	2-8°C	30 Sep 2021
Human α -Synuclein Calibrator	A0080167	$\leq -70^\circ\text{C}$	30 Sep 2020
Diluent 49	M0490004	$\leq -10^\circ\text{C}$	30 Sep 2019
Diluent 49	M0490004	$\leq -10^\circ\text{C}$	30 Sep 2019
Biotin Anti-hu α -Synuclein	D0081084	2-8°C	31 Mar 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in U-PLEX Plus	
20X MSD Wash Buffer	Not Kit Specific
Plate Seals	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>	01 MAR 2019