

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: K0081532

Expiration Date: 30 Apr 2022

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	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	α -Synuclein	8.00	80% - 120%	< 20%	6800	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	α -Synuclein	20X Calibrator	210000	pg/mL
		Human α -Synuclein Control 1	18400	pg/mL
		Human α -Synuclein Control 2	1740	pg/mL
		Human α -Synuclein Control 3	200	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:

α -Synuclein - Spot 1		
Conc. pg/mL	ECL Signal	%CV
10500	402,000	4.2
2630	79,900	4.0
656	16,400	4.3
164	3,760	4.3
41.0	1,050	4.8
10.3	324	2.8
2.56	145	4.3
0.00	85	4.9

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD Gold 96-Well Small Spot Streptavidin Plate	Z0021663	2-8°C	31 Oct 2022
Anti-hu α -Synuclein	D00U1065	$\leq -70^{\circ}\text{C}$	31 May 2022
Biotin Anti-hu α -Synuclein	B00U1065A	$\leq -70^{\circ}\text{C}$	30 Nov 2022
Human α -Synuclein Calibrator	A0080237	$\leq -70^{\circ}\text{C}$	30 Apr 2022
Diluent 49	M0490010	$\leq -10^{\circ}\text{C}$	30 Nov 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in U-PLEX Plus	
MSD Wash Buffer (20X), 100mL	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	Karen Hamilla	Quality		02 SEP 2020