

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: K0081330
Expiration Date: 31 Mar 2021

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	212000	pg/mL
		Human α -Synuclein Control 1	17000	pg/mL
		Human α -Synuclein Control 2	1610	pg/mL
		Human α -Synuclein Control 3	174	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
10600	261,000	2.4
2650	55,700	4.3
663	11,800	4.3
166	2,770	3.4
41.4	760	3.1
10.4	242	3.5
2.59	118	3.3
0.00	77	4.3

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD Gold 96-Well Small Spot Streptavidin Plate	Z0021482	2-8°C	30 Jun 2021
Anti-hu α -Synuclein	D00U0829	2-8°C	30 Sep 2021
Human α -Synuclein Calibrator	A0080211	\leq -70°C	30 Sep 2021
Diluent 49	M0490007	\leq -10°C	31 Oct 2021
Biotin Anti-hu α -Synuclein	B00U0830	2-8°C	31 Mar 2021
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		30 SEP 2019