

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: U-PLEX® and U-PLEX Plus Human α -Synuclein Kit

Kit Lot Number: K0081964

Expiration Date: 31 May 2023

Kit Catalog Numbers: K151WK-Series

Precision and Accuracy Specifications:

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

* See Additional Comments.

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

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	α -Synuclein	202000	pg/mL

Lot-Specific Components:


Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD™ 96-well Small Spot Streptavidin Plate	Z0021878	2-8°C	30 Jun 2024
Anti-hu α -Synuclein Antibody	D00U1872	\leq -70°C	31 Aug 2024
Biotin Anti-hu α -Synuclein Antibody	B00U1872	\leq -70°C	30 Sep 2024
Human α -Synuclein Calibrator	A0080272	\leq -70°C	31 May 2023
Diluent 49	M0490011	\leq -10°C	30 Nov 2023

Additional Comments:

	Assay	Spot	Specification	Comment
1	α -Synuclein	1	Control Conc.	% Control Recovery specification is 80-120%; the measured recovery for Control 1 & 2 was 123.2% & 120.6%

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 MAY 2022