

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: U-PLEX® and U-PLEX Plus Human α -Synuclein Kit
Kit Lot Number: K0081799
Expiration Date: 28 Feb 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%

		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 SECTOR Plate	Z0021706	2-8°C	28 Feb 2023
Anti-hu α -Synuclein	D00U1673	$\leq -70^\circ\text{C}$	31 Dec 2023
Human α -Synuclein Calibrator	A0080272	$\leq -70^\circ\text{C}$	31 May 2023
Diluent 49	M0490011	$\leq -10^\circ\text{C}$	30 Nov 2023
Biotin Anti-hu α -Synuclein	B00U1673	$\leq -70^\circ\text{C}$	31 Dec 2023

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		09 AUG 2021