



Meso Scale Diagnostics, LLC.

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Kit Lot Number: K0081716

Expiration Date: 30 Apr 2022

Kit Catalog Numbers: K151WK-Series

Precision and Accuracy 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 Calibrator:

Spot	Name	Conc.	Units
1	α -Synuclein	210000	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
Biotin Anti-hu α -Synuclein Antibody	B00U1498	$\leq -70^\circ\text{C}$	31 Aug 2023
Anti-hu α -Synuclein Antibody	D00U1306	$\leq -70^\circ\text{C}$	31 Jan 2023
Human α -Synuclein Calibrator	A0080237	$\leq -70^\circ\text{C}$	30 Apr 2022
Diluent 49	M0490010	$\leq -10^\circ\text{C}$	30 Nov 2022

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		06 MAY 2021