



CERTIFICATE OF ANALYSIS

Product Description: U-PLEX and U-PLEX Plus Human α -Synuclein Kit**Lot Number:** K0082832**Expiration Date:** 28 FEB 2027**Catalog Number:** K151WK-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD 96-Well Small Spot Streptavidin SECTOR Plate	Z0022570	2-8°C	29 FEB 2028
Diluent 49	M0490020	$\leq -10^{\circ}\text{C}$	28 FEB 2027
Anti-hu α -Synuclein	D00U2660	$\leq -70^{\circ}\text{C}$	28 FEB 2028
Biotin Anti-hu α -Synuclein	B00U2660	$\leq -70^{\circ}\text{C}$	31 JUL 2027
Human α -Synuclein Calibrator	A0080441	$\leq -70^{\circ}\text{C}$	29 FEB 2028

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

Precision, Accuracy, and Sensitivity 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	8.0	80% - 120%	< 20%	6800	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	α -Synuclein	214000	pg/mL

Dilution to top of Curve: 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 NOV 2025