

**CERTIFICATE OF ANALYSIS**

**Product Description:** U-PLEX® and U-PLEX Plus Human  $\alpha$ -Synuclein Kit  
**Lot Number:** K0082442  
**Expiration Date:** 31 JAN 2026  
**Catalog Number:** K151WK-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD™ 96-well Small Spot Streptavidin SECTOR Plate	Z0022202	2-8°C	31 JAN 2026
Diluent 49	M0490018	≤ -10°C	30 NOV 2026
Anti-hu $\alpha$ -Synuclein	D00U2482	≤ -70°C	31 DEC 2026
Biotin Anti-hu $\alpha$ -Synuclein	B00U2482	≤ -70°C	30 NOV 2026
Human $\alpha$ -Synuclein Calibrator	A0080399	≤ -70°C	30 NOV 2026

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

**Precision, Accuracy, and Sensitivity Specifications:**

Spot	Name	Calibrator		Human $\alpha$ -Synuclein Control 1		Human $\alpha$ -Synuclein Control 2		Human $\alpha$ -Synuclein Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	$\alpha$ -Synuclein	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

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

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	$\alpha$ -Synuclein	210000	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
<b>Review/Approval</b>	Karen Hamilla	Quality		23 JUL 2024