

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** U-PLEX® and U-PLEX Plus Human  $\alpha$ -Synuclein Kit  
**Kit Lot Number:** K0082011  
**Expiration Date:** 31 Dec 2023  
**Kit Catalog Numbers:** K151WK-Series

**Precision and Accuracy Specifications:**

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

		LLOQ			ULOQ		
Spot	Name	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	200000	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD™ 96-well Small Spot Streptavidin SECTOR Plate	Z0021925	2-8°C	30 Sep 2024
Biotin Anti-hu $\alpha$ -Synuclein	B00U1674	≤ -70°C	30 Sep 2024
Anti-hu $\alpha$ -Synuclein	D00U1673	≤ -70°C	31 Dec 2023
Human $\alpha$ -Synuclein Calibrator	A00U0270	≤ -70°C	30 Nov 2024
Diluent 49	M0490014	≤ -10°C	30 Sep 2024

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		08 JUL 2022