

**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** U-PLEX® and U-PLEX Plus Human  $\alpha$ -Synuclein Kit  
**Kit Lot Number:** K0082252  
**Expiration Date:** 31 May 2025  
**Kit Catalog Numbers:** K151WK-Series

**Precision and Accuracy 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	214000	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD™ 96-well Small Spot Streptavidin SECTOR Plate	Z0022045	2-8°C	31 May 2025
Anti-hu $\alpha$ -Synuclein	D00U2160	$\leq$ -70°C	31 Aug 2025
Human $\alpha$ -Synuclein Calibrator	A0080362	$\leq$ -70°C	31 Aug 2025
Diluent 49	M0490016	$\leq$ -10°C	30 Jun 2025
Biotin Anti-hu $\alpha$ -Synuclein	B00U2160	$\leq$ -70°C	31 Jul 2025

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		18 MAY 2023