

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** U-PLEX® and U-PLEX Plus Human  $\alpha$ -Synuclein Kit

**Kit Lot Number:** K0081903

**Expiration Date:** 31 May 2023

**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	202000	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Streptavidin Small Spot SECTOR Plate	Z0021837	2-8°C	29 Feb 2024
Anti-hu $\alpha$ -Synuclein	D00U1673	$\leq$ -70°C	31 Dec 2023
Human $\alpha$ -Synuclein Calibrator	A0080272	$\leq$ -70°C	31 May 2023
Diluent 49	M0490011	$\leq$ -10°C	30 Nov 2023
Biotin Anti-hu $\alpha$ -Synuclein	B00U1673	$\leq$ -70°C	31 Dec 2023

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		02 FEB 2022

