



Meso Scale Diagnostics, LLC.

### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** U-PLEX® and U-PLEX Plus Human  $\alpha$ -Synuclein Kit  
**Kit Lot Number:** K0081520  
**Expiration Date:** 30 Apr 2022  
**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	210000	pg/mL

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD 96-Well Small Spot Streptavidin Plate	Z0021663	2-8°C	31 Oct 2022
Anti-hu $\alpha$ -Synuclein	D00U1306	$\leq -70^\circ\text{C}$	31 Jan 2023
Human $\alpha$ -Synuclein Calibrator	A0080237	$\leq -70^\circ\text{C}$	30 Apr 2022
Diluent 49	M0490010	$\leq -10^\circ\text{C}$	30 Nov 2022
Biotin Anti-hu $\alpha$ -Synuclein	B00U1306	$\leq -70^\circ\text{C}$	31 May 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		20 JAN 2021