

MSD[®] α -Synuclein Calibrator



Ordering Information

MSD Customer Service
Phone: 1-240-314-2795
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com
www.mesoscale.com/support

Scientific Support

Phone: 1-240-314-2798
Email: ScientificSupport@mesoscale.com
www.mesoscale.com/support

Company Address

MESO SCALE DISCOVERY[®]
A division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA

www.mesoscale.com[®]

For Research Use Only.
Not for use in diagnostic
procedures.

Catalog #	C01WK-2
Contents	α -Synuclein Calibrator (1 vial; 30 μ L/vial)
Storage	$\leq -70^{\circ}\text{C}$

Summary and Intended Use

Human α -synuclein is a 140 amino acid protein abundantly expressed in the nervous system and genetically linked to Parkinson's disease. It is thought to maintain synaptic integrity and normal cellular homeostasis through synaptic vesicle recycling and neurotransmitter release modulation. The α -Synuclein Calibrator is intended for use as an assay standard for MSD assays. The utility of this product has not been evaluated in other applications.

Reagent

The α -Synuclein Calibrator is a full length recombinant human protein expressed in *E.coli* and is supplied as a 20X concentrated frozen liquid.

Storage and Handling

In order to maximize consistency in calibrator values across assay runs, the calibrator must be stored at the temperature recommended above. This calibrator can go through four freeze-thaw cycles. Unused material should be discarded after the fourth thaw.

The α -synuclein calibration curves are constructed based on the concentration of α -Synuclein Calibrator in each well. Refer to the assay-specific product insert for calibrator preparation instructions. Excess diluted material should be discarded after use.

Safety

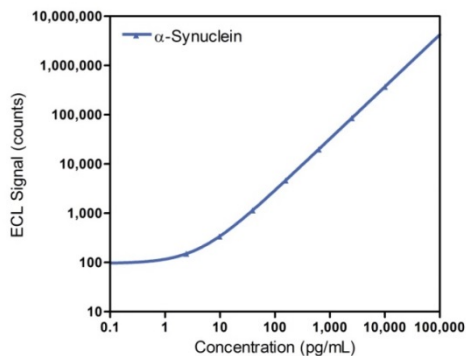
Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling calibrator. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines. Refer to the product-specific safety data sheet for additional safety and product handling information. The safety data sheet can be obtained from MSD Customer Service or at www.mesoscale.com.

Assignment of Values

The α -synuclein protein concentrations in a reference lot of α -Synuclein Calibrator were confirmed by spectrophotometry. For lot-specific calibrator concentrations, refer to the certificate of analysis (COA) supplied with the product.

Representative Data

The representative α -synuclein assay curve reported in this insert was determined using the U-PLEX[®] Human α -Synuclein Kit (Catalog #K151WKK-2). The reported data should be used only as a guide.



α -Synuclein		
Conc. (pg/mL)	Average Signal	%CV
10,000	370,227	4.6
2,500	85,449	4.3
625	19,778	3.9
156	4,634	3.0
39.1	1,142	2.5
9.77	337	3.2
2.44	151	3.3
0	95	6.7

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, www.mesoscale.com, SMALL SPOT (design), MSD, MSD (design), U-PLEX, and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. © 2016 Meso Scale Diagnostics, LLC. All rights reserved.

