

# MSD<sup>®</sup> Human $\alpha$ -Synuclein Control Pack



## Ordering Information

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

## Scientific Support

Phone: 1-240-314-2798  
Email: ScientificSupport@mesoscale.com  
[www.mesoscale.com/support](http://www.mesoscale.com/support)

## Company Address

MESO SCALE DISCOVERY<sup>®</sup>  
A division of  
Meso Scale Diagnostics, LLC.  
1601 Research Boulevard  
Rockville, MD 20850-3173 USA

[www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup>

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

<b>Catalog #</b>	C41WK-1
<b>Contents</b>	Human $\alpha$ -Synuclein Control 1 (5 vials*; 40 $\mu$ L/vial) Human $\alpha$ -Synuclein Control 2 (5 vials*; 40 $\mu$ L/vial) Human $\alpha$ -Synuclein Control 3 (5 vials*; 40 $\mu$ L/vial) *If supplied as part of a U-PLEX <sup>®</sup> Plus kit, quantity will match number of plates ordered.
<b>Storage</b>	$\leq -70^{\circ}\text{C}$

## Summary and Intended Use

The use of controls to monitor the analytical performance and reliability of test methods is an essential component of good laboratory practice.

The Human  $\alpha$ -Synuclein Control Pack consists of three levels of controls, each containing known concentrations of  $\alpha$ -synuclein protein. Human  $\alpha$ -Synuclein Controls 1, 2, and 3 are prepared by spiking recombinant calibrators into diluent. The controls are supplied as frozen liquid.

## Storage and Handling

To maximize the consistency of measured values across vials, the controls must be stored at the recommended temperature. Controls can go through four freeze-thaw cycles without significantly changing analyte levels. Discard unused material after the fourth thaw.

To use, dilute controls 8-fold in assay diluent. Add diluted control solutions to the MSD plate (coated with capture antibody) and assay as unknown samples. Discard excess diluted control material after use.

## Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling controls. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines. Additional product-specific safety information is available in the safety data sheet, which can be obtained from MSD Customer Service or at [www.mesoscale.com](http://www.mesoscale.com).

## Assignment of Control Values

The controls are provided to assess reproducibility of assay performance, and precision CVs are expected to be less than 20%. The certificate of analysis contains the concentrations of the controls measured at MSD, across three lots of the U-PLEX Human  $\alpha$ -Synuclein Kit (Catalog #K151WKK-1). Even with good laboratory practices, site-to-site differences may occur. Therefore, to set accuracy specifications, it is recommended that each lab establish its own nominal values and acceptance range for control concentrations.