

MSD[®] Neurodegeneration Control Pack 1 (6E10)



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 No.	C41LB-1
Contents	Neurodegeneration Control 1 (6E10) (5 vials,* 100 µL/vial) Neurodegeneration Control 2 (6E10) (5 vials,* 100 µL/vial) Neurodegeneration Control 3 (6E10) (5 vials,* 100 µL/vial) <small>*If supplied as part of a V-PLEX Plus kit, quantity will match number of plates ordered.</small>
Storage	≤-70°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 Neurodegeneration Control Pack 1 consists of three levels of controls, each containing known concentrations of Aβ38, Aβ40, and Aβ42 peptides and are intended for use as quality controls in MSD V-PLEX[®] assays that measure these analytes. Each vial contains enough materials for at least six measurements.

Reagent

Neurodegeneration Controls 1, 2, and 3 are prepared by spiking known levels of synthetic Aβ peptides into diluent. The controls are supplied as frozen solutions.

Storage and Handling

In order to maximize consistency of measurement across vials, the controls must be stored at the temperature recommended above. The controls can be subjected to three freeze–thaw cycles without significantly affecting the concentration of analytes. Unused material should be discarded after the third thaw.

For use as quality controls in MSD kits, we suggest handling Neurodegeneration Controls 1, 2, and 3 according to the recommended sample preparation method stated in the kit-specific product insert. Once prepared, the controls should be assayed within one hour. Excess diluted material should be discarded.

Assignment of Values

Neurodegeneration Control Pack 1 (6E10) analyte concentrations are determined using the following MSD kits:

V-PLEX Plus Kit Name	Kit Catalog Number
Aβ Peptide Panel 1 (6E10) Kit	K15200G
Aβ38 Peptide (6E10) Kit	K150SIG
Aβ40 Peptide (6E10) Kit	K150SKG
Aβ42 Peptide (6E10) Kit	K151LBG

Controls should be diluted according to the recommendations described in the respective product inserts. The controls are provided to assess reproducibility of assay performance. For the lot-specific concentration of each analyte in the control pack, refer to the certificate of analysis (COA) supplied with the kit or available at www.mesoscale.com. Even with good laboratory practices, site-to-site differences may occur. Therefore, to establish accuracy specifications, it is recommended that each lab should establish its own nominal values and acceptance range for the concentrations of the controls.

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 applicable safety data sheet(s) (SDS), which can be obtained from MSD Customer Service or at www.mesoscale.com.

MESO SCALE DISCOVERY, Meso Scale Diagnostics, MSD, V-PLEX, www.mesoscale.com, 96 WELL 4-SPOT (design), MSD (design), and Spot the Difference are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

The 6E10 antibody used in MSD Aβ assays is supplied by BioLegend (previously from Covance Research Products).

©2013, 2017, 2019, 2023 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-141-v5-2023oct