

MSD[®] Neurodegeneration A β Peptides

Calibrators



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[®]

Catalog #	Contents	Storage
C00NZ-2	A β 1-38 peptide (1 vial)	$\leq -70^{\circ}\text{C}$
C00OA-2	A β 1-40 peptide (1 vial)	$\leq -70^{\circ}\text{C}$
C01LB-2	A β 1-42 peptide (1 vial)	$\leq -70^{\circ}\text{C}$

Summary and Intended Use

MSD Neurodegeneration A β 1-38, A β 1-40, and A β 1-42 peptides are provided as individual solutions that may be blended for use in multiplex immunoassays. These peptide solutions were developed for use as calibrators for MSD A β assays. The utility of these products has not been evaluated in other applications.

Reagent

MSD's A β peptide calibrators are made from synthetic A β 1-38, A β 1-40, and A β 1-42 peptides and are supplied as frozen solutions. When used in MSD assays, the calibrators are diluted with assay diluent according to the recommendations in the assay product insert.

Storage and Handling

In order to maximize consistency in calibrator values across assay runs, the calibrator must be stored at the recommended temperature. When using these calibrators in MSD assays, please refer to the instructions included in the assay product insert. Once thawed, MSD recommends using within one hour. The calibrators can tolerate 3 freeze-thaw cycles. Excess diluted calibrator should be discarded.

Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling calibrators. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines. Refer to the safety data sheet, included in the product package and available from MSD Customer Service, for additional product safety and handling information.

Assignment of Values

A β peptide calibrator concentrations are calibrated against a reference calibrator generated at MSD. For lot-specific calibrator concentrations, refer to the certificate of analysis supplied with the calibrator.

Representative Data

The representative A β peptide calibration curves reported in this insert were determined using the MSD A β Peptide Panel 1 (6E10) Kit (catalog # K15200E). The reported data should be used only as a guide. The average calibrator signals provided below are specific to the lots tested.

A β 38		
Conc. (pg/mL)	Average Signal	%CV
10 000	648 926	5.0
2 500	100 970	5.5
625	9052	7.6
156	926	6.3
39.1	195	5.4
9.77	100	7.1
2.44	81	7.3
0	74	8.8

A β 40		
Conc. (pg/mL)	Average Signal	%CV
15 000	403 391	5.5
3 750	135 018	4.3
938	22 951	5.7
234	2590	6.1
58.6	410	6.5
14.6	134	5.1
3.66	88	8.2
0	74	7.0

A β 42		
Conc. (pg/mL)	Average Signal	%CV
1 450	725 654	3.3
363	182 208	3.2
91	31 037	3.4
23	4551	4.2
5.7	917	4.9
1.4	268	4.6
0.35	134	6.1
0	90	6.8

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

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, www.mesoscale.com, 96 WELL 1, 4, 7, & 10-SPOT (designs), MSD (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. © 2013 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-187-v1-2013Aug