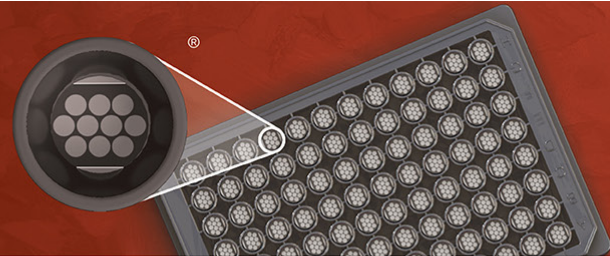


MSD® Metabolic Panel 1 Calibrator Blend



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: 1-240-314-2795
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 #	C0325
Contents	Metabolic Panel 1 Calibrator Blend (1 vial; lyophilized)
Storage	2–8 °C

Summary and Intended Use

The Metabolic Panel 1 Calibrator Blend is a lyophilized blend of seven metabolic proteins: insulin, C-Peptide, GLP-1, GIP, glucagon, PP, and leptin. The Metabolic Panel 1 Calibrator Blend was developed for use as an assay calibrator for MSD assays in the V-PLEX® Plus and V-PLEX Metabolic Panel 1 (human) Kits (catalog # K15325D, K15325G); V-PLEX Plus and V-PLEX Metabolic Panel 1 (NHP) Kits (catalog # K15332D, K15332G); V-PLEX Plus and V-PLEX Metabolic Panel 1 (mouse) Kits (catalog # K15333D, K15333G); V-PLEX Plus and V-PLEX Metabolic Panel 1 (rat) Kits (catalog # K15334D, K15334G); and V-PLEX Plus and V-PLEX Metabolic Panel 1 (canine) Kits (catalog # K15335D, K15335G). The utility of this product has not been evaluated in other kits beside the one mentioned above.

Reagent

Metabolic Panel 1 Calibrator Blend includes full-length *E. coli* recombinant proteins (insulin, glucagon, and leptin); and synthetic peptides (C-Peptide, GIP (1-42), PP, and amidated GLP-1 (7-36)). The Metabolic Panel 1 Calibrator Blend is supplied as a lyophilized powder that yields top-of-curve calibrator concentrations upon reconstitution in 1 mL of assay diluent (Diluent 56; catalog # R50BR).

Storage and Handling

In order to maximize consistency in calibrator values across assay runs, the calibrator must be stored at the recommended temperature and handled per the instructions provided in this document. Once reconstituted in 1 mL of assay diluent (Diluent 56; catalog # R50BR); the calibrator is stable at $\leq -70^{\circ}\text{C}$ and can tolerate three freeze-thaw cycles.

Metabolic Panel 1 calibration curves are constructed based on the concentration of calibrator in each well. Refer to the assay-specific product insert for calibrator preparation instructions. Excess diluted calibrator should be discarded after use.

For the lot-specific concentration of the calibrator, refer to the certificate of analysis (COA) supplied with the kit or available at www.mesoscale.com.

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. Additional product-specific safety information is available in the safety data sheet (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 SMALL-SPOT and 10-SPOT (designs), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

©2020 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-914-v1-2020Mar