

MSD[®] GLP-1 (active) 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 #	C0266-1
Contents	GLP-1 (active) Calibrator (1 vial; lyophilized)
Storage	2-8°C

Summary and Intended Use

The GLP-1 (active) Calibrator was developed for use as an assay calibrator for MSD assays in the V-PLEX[®] GLP-1 Active and V-PLEX GLP-1 Active Plus Kits (catalog # K1503OD, K1503OG). The utility of this product has not been evaluated in other applications.

Reagent

GLP-1 (active) Calibrator includes synthetic amidated GLP-1 (7-36) peptide. The GLP-1 (active) Calibrator is supplied as a lyophilized powder that yields top-of-curve calibrator concentration upon reconstitution in 1 mL of Metabolic Assay Working Solution (see the V-PLEX GLP-1 Active Kit product insert for details).

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 Metabolic Assay Working Solution, the calibrator is stable at $\leq -70^{\circ}\text{C}$ and can tolerate three freeze–thaw cycles.

GLP-1 (active) assay 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, MSD (design), SMALL SPOT (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. ©2018 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-629-v1-2018Jan