

Angiogenesis Control Pack 1



www.mesoscale.com®

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: 1-301-947-2025
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

Catalog No.	C4666-1
Contents	Angiogenesis Control 1 (5 vials; * 150 µL/vial) Angiogenesis Control 2 (5 vials; * 150 µL/vial) Angiogenesis Control 3 (5 vials; * 150 µL/vial) *Supplied as part of a V-PLEX® Plus kit, quantity will match number of plates ordered.
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 Angiogenesis Control Pack 1 consists of 3 levels of controls, each containing known concentrations of all human angiogenesis proteins that are detected by the V-PLEX and V-PLEX Plus Angiogenesis Panel 1 (human) Kits (catalog No. K15190D and K15190G). The controls span the linear ranges of the assays in the V-PLEX and V-PLEX Plus Angiogenesis Panel 1 (human) Kits and are designed to test the analytical range of the assays. Angiogenesis Controls 1, 2, and 3 are prepared by spiking recombinant calibrators into a nonmatrix-based diluent. The controls are supplied frozen.

Note: The control pack (catalog No. C4190-1) was discontinued and replaced by a control pack containing new recombinant VEGF-C protein (catalog No. C4666-1).

Storage and Handling

To maximize the consistency of measured values across vials, the controls must be stored at the temperature recommended above. The controls can tolerate 3 freeze-thaw cycles. Discard unused material after the third freeze-thaw cycle.

To use, thaw each vial of controls on wet ice for a minimum of 30 minutes. Dilute controls 2-fold in assay diluent (Diluent 7). Add diluted control solution directly to the MSD Angiogenesis Panel 1 (human) plate and assay as unknown samples. Discard unused diluted control material.

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.

Assignment of Control Values

The controls are provided to assess reproducibility of assay performance, and precision CVs are expected to be less than 25%. The certificate of analysis contains the concentrations of the controls measured at MSD across three lots. 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 control concentrations.

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

MESO SCALE DISCOVERY, Meso Scale Diagnostics, MSD, www.mesoscale.com, MSD (design), V-PLEX, 96 WELL 10-SPOT (design), and Spot the Difference are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
©2014, 2021, 2023 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-1169-v1-2023Sep