

# Angiogenesis Panel 1 (human) Calibrator Blend



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](http://www.mesoscale.com/support)

## Scientific Support

Phone: 1-301-947-2025  
Email: ScientificSupport@mesoscale.com  
[www.mesoscale.com/support](http://www.mesoscale.com/support)

## Company Address

MESO SCALE DISCOVERY®  
A division of  
Meso Scale Diagnostics, LLC.  
1601 Research Boulevard  
Rockville, MD 20850-3173  
USA

<b>Catalog No.</b>	C0666-2
<b>Contents</b>	Angiogenesis Panel I (human) Calibrator Blend (1 vial; 60 µL)
<b>Storage</b>	≤-70°C

## Summary and Intended Use

The Angiogenesis Panel 1 (human) Calibrator Blend is a frozen, liquid combination of 7 recombinant human proteins: VEGF-A, VEGF-C, VEGF-D, Tie-2, Flt-1, PlGF, and FGF (basic). Angiogenesis Panel 1 (human) Calibrator Blend was developed for use as an assay calibrator for MSD assays in the V-PLEX® Angiogenesis Panel 1 (human) Kit (catalog No. K15190D-1). The utility of this product has not been evaluated in other applications.

**Note:** The calibrator blend (catalog No. C0190-2) was discontinued and replaced by a calibrator blend containing new recombinant VEGF-C protein (catalog No. C0666-2).

## Reagent

Angiogenesis Panel 1 (human) Calibrator Blend includes full-length recombinant proteins for each analyte. Angiogenesis Panel 1 (human) calibrator blend is supplied as a concentrated liquid that yields top-of-curve calibrator concentration upon 20-fold dilution in assay diluent (Diluent 7, catalog No. R54BB).

## 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. The calibrator blend is stable at ≤-70 °C and can tolerate 1 freeze-thaw cycle.

Angiogenesis Panel 1 (human) 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.

## 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.

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, 2023 Meso Scale Diagnostics, LLC. All rights reserved.



MK-DS-1170-v1-2023Sep