

**CERTIFICATE OF ANALYSIS**


Product Description: Angiogenesis Panel 1 (human) Calibrator Blend
Catalog Number: C0666-2
Lot Number: A0080415
Expiration Date: 30 SEP 2027
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
VEGF-A	39400	pg/mL
VEGF-C	686000	pg/mL
VEGF-D	582000	pg/mL
Tie-2	1864000	pg/mL
VEGFR-1/Flt-1	204000	pg/mL
PlGF	24400	pg/mL
FGF (basic)	51200	pg/mL

Dilution to Top of Curve: 20 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		22 MAY 2025