



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Angiogenesis Panel 1 (human) Calibrator Blend

**Catalog Number:** C0190-2

**Lot Number:** A0080260

**Expiration Date:** 31 DEC 2022

**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
VEGF	40200	pg/mL
VEGF-C	536000	pg/mL
VEGF-D	612000	pg/mL
Tie-2	2020000	pg/mL
Flt-1	204000	pg/mL
PlGF	24000	pg/mL
bFGF	54400	pg/mL

**Dilution to Top of Curve:** 20X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		22 JUN 2020