



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Angiogenesis Panel 1 (human) Calibrator Blend

Catalog Number: C0190-2

Lot Number: A0080213

Expiration Date: 31 MAY 2022

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
VEGF-A	46700	pg/mL
VEGF-C	534000	pg/mL
VEGF-D	551000	pg/mL
Tie-2	1920000	pg/mL
Flt-1	187000	pg/mL
PlGF	22500	pg/mL
bFGF	50200	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
Review/Approval	Karen Hamilla	Quality		19 DEC 2019