



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Angiogenesis Panel 1 (human) Calibrator Blend

Catalog Number: C0190-2

Lot Number: A0080185

Expiration Date: 31 MAR 2021

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

Concentration of Materials Tested:

	Assay						
	hu VEGF	hu VEGF-C	hu VEGF-D	hu Tie-2	hu Flt-1	hu PIGF	hu bFGF
Calibrator (pg/mL)	49 000	492 000	568 000	1 840 000	196 600	24 200	49 800

Concentration: 20X

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	29 NOV 2018