



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Angiogenesis Panel 1 (human) Calibrator Blend

Catalog Number: C0190-2

Lot Number: A0080104

Expiration Date: 31 MAR 2019

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 PlGF	hu bFGF
Reconstituted Calibrator (pg/mL)	43000	414000	476000	1414000	162800	24000	37000

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 JUN 2017