



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Angiogenesis Control Pack 1

**Catalog Number:** C4190-1

**Lot Number:** CP000077

**Expiration Date:** 30 NOV 2020

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

### Concentration of Materials Tested:

Units = pg/mL	Assay						
	hu VEGF	hu VEGF-C	hu VEGF-D	hu Tie-2	hu Flt-1	hu PIGF	hu bFGF
Control 1 (A00C0594)	444	5710	6720	20 000	2310	260	559
Control 2 (A00C0595)	89.4	1560	896	6480	291	61.6	62.7
Control 3 (A00C0596)	16.6	380	92.2	444	27.0	12.2	5.88

**Concentration:** 2x

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Control 1	A00C0594	$\leq -70^{\circ}\text{C}$	30 NOV 2020
Angiogenesis Control 2	A00C0595	$\leq -70^{\circ}\text{C}$	30 NOV 2020
Angiogenesis Control 3	A00C0596	$\leq -70^{\circ}\text{C}$	30 NOV 2020

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		09 APR 2019