



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Control Pack Description:** Angiogenesis Control Pack 1

**Catalog Number:** C4190-1

**Lot Number:** CP000098

**Expiration Date:** 31 MAY 2022

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
VEGF	538	99.8	24.8	pg/mL
VEGF-C	6950	1840	614	pg/mL
VEGF-D	6530	1060	133	pg/mL
Tie-2	20200	6770	763	pg/mL
Flt-1	1800	214	32.0	pg/mL
PlGF	231	53.5	15.9	pg/mL
bFGF	567	64.7	7.85	pg/mL

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Control 1	A00C0667	≤ -70°C	31 MAY 2022
Angiogenesis Control 2	A00C0668	≤ -70°C	31 MAY 2022
Angiogenesis Control 3	A00C0669	≤ -70°C	31 MAY 2022

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		30 JUL 2020