



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C4190-1

**Lot Number:** CP000108

**Expiration Date:** 31 MAR 2023


**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
VEGF	461	99.2	20.7	pg/mL
VEGF-C	5220	1770	510	pg/mL
VEGF-D	5980	1130	144	pg/mL
Tie-2	16200	7590	1040	pg/mL
Flt-1	2250	253	35.0	pg/mL
PlGF	232	52.5	18.1	pg/mL
bFGF	529	68.3	8.73	pg/mL

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Control 1	A00C0720	≤ -70°C	31 MAR 2023
Angiogenesis Control 2	A00C0741	≤ -70°C	31 MAR 2023
Angiogenesis Control 3	A00C0742	≤ -70°C	31 MAR 2023

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Carissa B. Hubbard	Quality		26 OCT 2020