



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Angiogenesis Control Pack 1

Catalog Number: C4190-1

Lot Number: CP000117

Expiration Date: 30 SEP 2023

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
VEGF	547	113	21.8	pg/mL
VEGF-C	6070	1780	539	pg/mL
VEGF-D	8020	1400	245	pg/mL
Tie-2	26600	6820	1640	pg/mL
Flt-1	2320	290	35.6	pg/mL
PlGF	284	60.1	13.8	pg/mL
bFGF	602	69.9	8.13	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Control 1	A00C0779	$\leq -70^{\circ}\text{C}$	30 SEP 2023
Angiogenesis Control 2	A00C0780	$\leq -70^{\circ}\text{C}$	30 SEP 2023
Angiogenesis Control 3	A00C0781	$\leq -70^{\circ}\text{C}$	30 SEP 2023

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 JUN 2021