



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Angiogenesis Control Pack 1

Catalog Number: C4190-1

Lot Number: CP000122

Expiration Date: 31 DEC 2023

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
VEGF-A	605	110	18.1	pg/mL
VEGF-C	5930	1370	413	pg/mL
VEGF-D	8730	1580	282	pg/mL
Tie-2	32600	7960	1820	pg/mL
Flt-1	2220	131	36.9	pg/mL
PIGF	290	70.3	16.2	pg/mL
bFGF	610	78.7	7.88	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Control 1	A00C0793	≤ -70°C	31 DEC 2023
Angiogenesis Control 2	A00C0794	≤ -70°C	31 DEC 2023
Angiogenesis Control 3	A00C0795	≤ -70°C	31 DEC 2023

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 SEP 2021