



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Angiogenesis Control Pack 1

Catalog Number: C4190-1

Lot Number: CP000113

Expiration Date: 31 MAR 2023

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
VEGF	556	114	20.7	pg/mL
VEGF-C	7070	1930	510	pg/mL
VEGF-D	8300	1460	144	pg/mL
Tie-2	28400	7450	1040	pg/mL
Flt-1	2420	297	35.0	pg/mL
PlGF	259	61.2	18.1	pg/mL
bFGF	545	60.3	8.73	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Control 1	A00C0768	≤ -70°C	31 JUL 2023
Angiogenesis Control 2	A00C0769	≤ -70°C	31 JUL 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
Review/Approval	Karen Hamilla	Quality		09 MAR 2021