

**Extension Memo**

Product Description: Angiogenesis Control Pack 1
Lot Number: CP000146
New Expiration Date: 30 SEP 2027
Catalog Number(s): C4666-1

Based on additional testing, the original date of expiration for this control pack has been extended by 24 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Angiogenesis Control 1	A00C0886	30 SEP 2025	30 SEP 2027
Angiogenesis Control 2	A00C0887	30 SEP 2025	30 SEP 2027
Angiogenesis Control 3	A00C0888	30 SEP 2025	30 SEP 2027

Function	Name	Signature	Date
Quality	Karen Hamilla		28 Jul 2026



CERTIFICATE OF ANALYSIS

Control Pack Description: Angiogenesis Control Pack 1
Catalog Number: C4666-1
Lot Number: CP000146
Expiration Date: 30 SEP 2025

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
VEGF-A	526	147	44	pg/mL
VEGF-C	6830	2600	1090	pg/mL
VEGF-D	8780	2110	547	pg/mL
Tie-2	31300	11100	3960	pg/mL
VEGFR-1/Flt-1	2660	471	77.6	pg/mL
PlGF	293	91.1	32.5	pg/mL
FGF (basic)	586	95.0	16.8	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Control 1	A00C0886	$\leq -70^{\circ}\text{C}$	30 SEP 2025
Angiogenesis Control 2	A00C0887	$\leq -70^{\circ}\text{C}$	30 SEP 2025
Angiogenesis Control 3	A00C0888	$\leq -70^{\circ}\text{C}$	30 SEP 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 MAY 2023