



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Angiogenesis Control Pack

**Catalog Number:** C4190-1

**Lot Number:** CP000049

**Expiration Date:** 31 JAN 2019

**Concentration of Materials Tested:**

Units = pg/mL	Assay						
	hu VEGF	hu VEGF-C	hu VEGF-D	hu Tie-2	hu Flt-1	hu PlGF	hu bFGF
Angiogenesis Control 1 (A00C0464)	507	5520	5090	15790	2330	239	544
Angiogenesis Control 2 (A00C0465)	108	1550	1210	7520	274	48.6	62.6
Angiogenesis Control 3 (A00C0423)	20.6	457	234	1380	29.9	15.8	7.37

Control concentrations were assigned across a minimum of three kit lots.

Angiogenesis Control Pack 1 Components			
Description	Lot Number	Storage Temperature	Expiration Date
Angiogenesis Control 1	A00C0464	≤ -70°C	31 JUL 2019
Angiogenesis Control 2	A00C0465	≤ -70°C	31 JUL 2019
Angiogenesis Control 3	A00C0423	≤ -70°C	31 JAN 2019

**Statement:** The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	03 APR 2017