



Human Angiogenesis ELISA Strip II Protein Standards

Catalog Number EA-1042

(For Research Use Only)

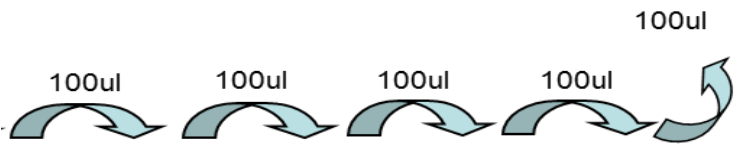
<u>Tube</u>	<u>Protein</u>	<u>Stock Conc.</u>	<u>Volume</u>
#1	Human PDGF-BB	400ng/ml	10ul
#2	Human PIGF-1	400ng/ml	10ul
#3	Human β -NGF	400ng/ml	10ul
#4	Human SCF	400ng/ml	10ul
#5	Human MCP-1	400ng/ml	10ul
#6	Human MIP-1 α	400ng/ml	10ul
#7	Human IL-2	400ng/ml	10ul
#8	Human IL-4	400ng/ml	10ul

Preparation of protein standard dilutions

1. Add 200ul of Diluent buffer the wells of the first strip, and add 100ul of Diluent buffer to the wells of the rest of the strips according to the following table.
2. Add appropriate amount of protein standards to the first strip according to the table 1.
3. Use multi-channel pipette to mix the dilutions and transfer 100ul to the next dilution wells.
4. Repeat the transfer until 5th strip. Trash 100ul from 5th strip after mixing. Do not transfer to the 6th Strip

Note: Substrate incubation time may vary due to different antibodies reactivity. Stronger signals (Strong blue color) could be stopped early after 5 minutes. Weaker signals should be incubated for 10-30 minutes. Always stop protein standards along with samples from the same row at the same time.

Table 1: Dilution of protein standards



<u>Protein</u>	<u>Added to 1st Strip</u>	<u>1st Strip</u>	<u>2nd Strip</u>	<u>3rd Strip</u>	<u>4th Strip</u>	<u>5th Strip</u>	<u>6th Strip</u>
Human PDGF-BB	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human PIGF-1	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human β -NGF	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human SCF	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human MCP-1	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human MIP-1 α	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human IL-2	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human IL-4	4ul	200ul	100ul	100ul	100ul	100ul	100ul
			1:2	1:4	1:8	1:16	Blank

Table 2: Concentrations of Protein Standard Dilutions

<u>Protein</u>		<u>1:2</u>	<u>1:4</u>	<u>1:8</u>	<u>1:16</u>	<u>Blank</u>
Human PDGF-BB	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human PIGF-1	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human β -NGF	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human SCF	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human MCP-1	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human MIP-1 α	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human IL-2	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human IL-4	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank