



Human Inflammation ELISA Strip Protein Standards

Catalog Number EA-1032

(For Research Use Only)

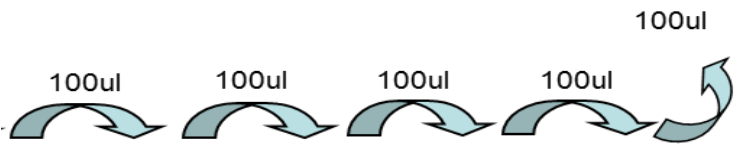
<u>Tube</u>	<u>Protein</u>	<u>Stock Conc.</u>	<u>Volume</u>
#1	Human TNF α	400ng/ml	10ul
#2	Human IFN γ	400ng/ml	10ul
#3	Human G-CSF	400ng/ml	10ul
#4	Human GM-CSF	400ng/ml	10ul
#5	Human IL-1 α	400ng/ml	10ul
#6	Human IL-8	400ng/ml	10ul
#7	Human IP-10	400ng/ml	10ul
#8	Human Rantes	400ng/ml	10ul

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>
		200ul	100ul	100ul	100ul	100ul	100ul
Human TNF α	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human IFN γ	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human G-CSF	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human GM-CSF	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human IL-1 α	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human IL-8	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human IP-10	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Human Rantes	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 TNF α	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human IFN γ	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human G-CSF	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human GM-CSF	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human IL-1 α	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human IL-8	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human IP-10	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Human Rantes	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank