



## Mouse Inflammation ELISA Strip Protein Standards

Catalog Number EA-1052

(For Research Use Only)

<u>Tube</u>	<u>Protein</u>	<u>Stock Conc.</u>	<u>Volume</u>
#1	Mouse IL-1 $\alpha$	400ng/ml	10ul
#2	Mouse IL-1 $\beta$	400ng/ml	10ul
#3	Mouse G-CSF	400ng/ml	10ul
#4	Mouse GM-CSF	400ng/ml	10ul
#5	Mouse MCP-1	400ng/ml	10ul
#6	Mouse MIP-1 $\alpha$	400ng/ml	10ul
#7	Mouse SCF	400ng/ml	10ul
#8	Mouse 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 5<sup>th</sup> strip. Trash 100ul from 5<sup>th</sup> strip after mixing. Do not transfer to the 6<sup>th</sup> 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>
Mouse IL-1 $\alpha$	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Mouse IL-1 $\beta$	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Mouse G-CSF	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Mouse GM-CSF	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Mouse MCP-1	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Mouse MIP-1 $\alpha$	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Mouse SCF	4ul	200ul	100ul	100ul	100ul	100ul	100ul
Mouse Rantes	4ul	200ul	100ul	100ul	100ul	100ul	100ul
			<b>1:2</b>	<b>1:4</b>	<b>1:8</b>	<b>1:16</b>	<b>Blank</b>

**Table 2: Concentrations of Protein Standard Dilutions**

<b><u>Protein</u></b>		<b><u>1:2</u></b>	<b><u>1:4</u></b>	<b><u>1:8</u></b>	<b><u>1:16</u></b>	<b><u>Blank</u></b>
Mouse IL-1 $\alpha$	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Mouse IL-1 $\beta$	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Mouse G-CSF	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Mouse GM-CSF	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Mouse MCP-1	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Mouse MIP-1 $\alpha$	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Mouse SCF	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank
Mouse Rantes	8ng/ml	4ng/ml	2ng/ml	1ng/ml	0.5ng/ml	Blank