

## JBScreen Membrane 2

Cat.-No: CS-302L

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
<b>A 1</b>	12% w/v PEG 2000 MME	None	50 mM Tris-HCl	7.5	500 mM Sodium Chloride
<b>A 2</b>	10% w/v PEG 3350	None	50 mM Sodium Citrate	5.6	150 mM Sodium Chloride
<b>A 3</b>	2% w/v PEG 4000	None	50 mM Tris-HCl	7.5	None
<b>A 4</b>	5% w/v PEG 4000	None	None		None
<b>A 5</b>	5% w/v PEG 4000	None	None		100 mM Potassium Chloride
<b>A 6</b>	5% w/v PEG 4000	10% w/v Glycerol	50 mM MES Sodium Salt	6.5	100 mM Sodium Chloride
<b>B 1</b>	5% w/v PEG 4000	None	50 mM Sodium Phosphate	6.7	None
<b>B 2</b>	10% w/v PEG 4000	None	50 mM Tris-HCl	8.5	500 mM Sodium Chloride
<b>B 3</b>	12% w/v PEG 4000	None	100 mM ADA	6.5	100 mM Lithium Sulfate
<b>B 4</b>	12% w/v PEG 4000	None	50 mM Sodium Phosphate	6.8	None
<b>B 5</b>	12% w/v PEG 4000	20% w/v Glycerol	50 mM MOPS	7.0	500 mM Potassium Chloride
<b>B 6</b>	15% w/v PEG 4000	None	10 mM Tris-HCl	7.5	100 mM Lithium Chloride
<b>C 1</b>	20% w/v PEG 4000	None	100 mM Bis-Tris	7.0	500 mM Sodium Chloride
<b>C 2</b>	20% w/v PEG 4000	None	100 mM Sodium Phosphate	7.0	500 mM Sodium Chloride
<b>C 3</b>	20% w/v PEG 4000	150 mM Zinc Acetate	50 mM Tris-HCl	7.5	50 mM Zinc Chloride
<b>C 4</b>	22% w/v PEG 4000	None	50 mM Tricine	8.0	None
<b>C 5</b>	22% w/v PEG 4000	None	50 mM Tris-HCl	8.5	500 mM Sodium Chloride
<b>C 6</b>	30% w/v PEG 4000	None	None		None
<b>D 1</b>	10% w/v PEG 5000 MME	None	100 mM Sodium Citrate	5.6	100 mM Magnesium Acetate
<b>D 2</b>	5% w/v PEG 6000	None	None		100 mM Magnesium Sulfate
<b>D 3</b>	10% w/v PEG 6000	150 mM Zinc Acetate	50 mM Tris-HCl	7.5	50 mM Zinc Chloride
<b>D 4</b>	15% w/v PEG 6000	None	50 mM Sodium Succinate	6.5	None
<b>D 5</b>	12% w/v PEG 8000	10% w/v MPD	None		25 mM Potassium dihydrogen Phosphate
<b>D 6</b>	8% w/v PEG 10000	None	100 mM Sodium Citrate	5.6	100 mM Magnesium Acetate