

## JBScreen Classic 2

Cat.-No: CS-102L

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
<b>A 1</b>	4 % w/v PEG 4000	None	100 mM Sodium Acetate	4.6	None
<b>A 2</b>	8 % w/v PEG 4000	None	None		None
<b>A 3</b>	8 % w/v PEG 4000	None	100 mM Sodium Acetate	4.6	None
<b>A 4</b>	10 % w/v PEG 4000	None	100 mM MES Sodium Salt	6.5	200 mM Magnesium Chloride
<b>A 5</b>	12 % w/v PEG 4000	None	100 mM HEPES Sodium Salt	7.5	100 mM Sodium Acetate
<b>A 6</b>	12 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	None
<b>B 1</b>	16 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate
<b>B 2</b>	16 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	200 mM Sodium Acetate
<b>B 3</b>	16 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	200 mM Magnesium Chloride
<b>B 4</b>	18 % w/v PEG 4000	None	100 mM Sodium Acetate	4.6	None
<b>B 5</b>	20 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate
<b>B 6</b>	20 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	200 mM Calcium Chloride
<b>C 1</b>	22 % w/v PEG 4000	None	100 mM HEPES Sodium Salt	7.5	100 mM Sodium Acetate
<b>C 2</b>	25 % w/v PEG 4000	None	100 mM Sodium Acetate	4.6	None
<b>C 3</b>	25 % w/v PEG 4000	None	100 mM MES Sodium Salt	6.5	200 mM Magnesium Chloride
<b>C 4</b>	25 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	200 mM Calcium Chloride
<b>C 5</b>	30 % w/v PEG 4000	None	None		None
<b>C 6</b>	30 % w/v PEG 4000	None	100 mM Sodium Acetate	4.6	100 mM Magnesium Chloride
<b>D 1</b>	30 % w/v PEG 4000	None	100 mM MES Sodium Salt	6.5	None
<b>D 2</b>	30 % w/v PEG 4000	None	100 mM HEPES Sodium Salt	7.5	200 mM Calcium Chloride
<b>D 3</b>	30 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate
<b>D 4</b>	30 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	200 mM Sodium Acetate
<b>D 5</b>	30 % w/v PEG 4000	None	100 mM Tris-HCl	8.5	200 mM Magnesium Chloride
<b>D 6</b>	35 % w/v PEG 4000	None	None		None