

## JBScreen Classic 1

Cat.-No: CS-101L

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
<b>A 1</b>	15 % w/v PEG 400	None	100 mM Sodium Acetate	4.6	100 mM Calcium Chloride
<b>A 2</b>	15 % w/v PEG 400	None	100 mM MES Sodium Salt	6.5	None
<b>A 3</b>	15 % w/v PEG 400	None	100 mM HEPES Sodium Salt	7.5	200 mM Magnesium Chloride
<b>A 4</b>	15 % w/v PEG 400	None	100 mM Tris-HCl	8.5	200 mM Sodium Citrate
<b>A 5</b>	25 % w/v PEG 400	None	100 mM Sodium Acetate	4.6	100 mM Magnesium Chloride
<b>A 6</b>	25 % w/v PEG 400	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate
<b>B 1</b>	28 % w/v PEG 400	None	100 mM HEPES Sodium Salt	7.5	200 mM Calcium Chloride
<b>B 2</b>	30 % w/v PEG 400	None	100 mM Sodium Acetate	4.6	100 mM Calcium Chloride
<b>B 3</b>	30 % w/v PEG 400	None	100 mM MES Sodium Salt	6.5	100 mM Sodium Acetate
<b>B 4</b>	30 % w/v PEG 400	None	100 mM MES Sodium Salt	6.5	100 mM Magnesium Chloride
<b>B 5</b>	30 % w/v PEG 400	None	100 mM HEPES Sodium Salt	7.5	200 mM Magnesium Chloride
<b>B 6</b>	30 % w/v PEG 400	None	100 mM Tris-HCl	8.5	200 mM Sodium Citrate
<b>C 1</b>	30 % w/v PEG 550 MME	None	100 mM Bicine	9.0	100 mM Sodium Chloride
<b>C 2</b>	25 % w/v PEG 550 MME	None	100 mM MES Sodium Salt	6.5	10 mM Zinc Sulfate
<b>C 3</b>	25 % w/v PEG 1000	None	100 mM HEPES Sodium Salt	7.5	None
<b>C 4</b>	30 % w/v PEG 1000	None	100 mM Tris-HCl	8.5	None
<b>C 5</b>	15 % w/v PEG 1500	None	None		None
<b>C 6</b>	20 % w/v PEG 1500	None	100 mM HEPES Sodium Salt	7.5	None
<b>D 1</b>	30 % w/v PEG 1500	None	None		None
<b>D 2</b>	20 % w/v PEG 2000 MME	None	100 mM Tris-HCl	8.5	10 mM Nickel(II) Chloride
<b>D 3</b>	25 % w/v PEG 2000 MME	None	None		None
<b>D 4</b>	30 % w/v PEG 2000 MME	None	100 mM MES Sodium Salt	6.5	100 mM Sodium Acetate
<b>D 5</b>	20 % w/v PEG 3000	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Acetate
<b>D 6</b>	30 % w/v PEG 3000	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate