

## JBScreen JCSG++ 1

Cat.-No: CS-151

Number	Precipitant 1	Precipitant 2	Buffer	pH	Salt	pH
<b>A 1</b>	50% v/v PEG 400	None	100 mM Sodium Acetate	4.5	200 mM Lithium Sulfate	
<b>A 2</b>	20% w/v PEG 3000	None	100 mM Sodium Citrate	5.5	None	
<b>A 3</b>	20% w/v PEG 3350	None	None		200 mM di-Ammonium hydrogen Citrate	5.0
<b>A 4</b>	30% v/v MPD	None	100 mM Sodium Acetate	4.6	20 mM Calcium Chloride	
<b>A 5</b>	20% w/v PEG 3350	None	None		200 mM Magnesium Formate	5.9
<b>A 6</b>	20% w/v PEG 1000	None	100 mM Phosphate / Citrate	4.2	200 mM Lithium Sulfate	
<b>B 1</b>	20% w/v PEG 8000	None	100 mM CHES	9.5	None	
<b>B 2</b>	20% w/v PEG 3350	None	None		200 mM Ammonium Formate	6.6
<b>B 3</b>	20% w/v PEG 3350	None	None		200 mM Ammonium Chloride	6.3
<b>B 4</b>	20% w/v PEG 3350	None	None		200 mM Potassium Formate	7.3
<b>B 5</b>	50% v/v MPD	None	100 mM Tris-HCl	8.5	200 mM Ammonium dihydrogen Phosphate	
<b>B 6</b>	20% w/v PEG 3350	None	None		200 mM Potassium Nitrate	6.9
<b>C 1</b>	None	None	100 mM Citric Acid	4.0	800 mM Ammonium Sulfate	
<b>C 2</b>	20% w/v PEG 3550	None	None		200 mM Sodium Thiocyanate	6.9
<b>C 3</b>	20% w/v PEG 6000	None	100 mM Bicine	9.0	None	
<b>C 4</b>	10% w/v PEG 8000	8% v/v Ethylene Glycol	100 mM HEPES Sodium Salt	7.5	None	
<b>C 5</b>	40% v/v MPD	5% w/v PEG 8000	100 mM MES Sodium Salt	6.5	None	
<b>C 6</b>	40% v/v Ethanol	5% w/v PEG 1000	100 mM Phosphate / Citrate	4.2	None	
<b>D 1</b>	8% w/v PEG 4000	None	100 mM Sodium Acetate	4.6	None	
<b>D 2</b>	10% w/v PEG 8000	None	100 mM Tris-HCl	7.0	200 mM Magnesium Chloride	
<b>D 3</b>	20% w/v PEG 6000	None	100 mM Citric Acid	5.0	None	
<b>D 4</b>	50% v/v PEG 200	None	100 mM MES Sodium Salt	6.5	200 mM Magnesium Chloride	
<b>D 5</b>	None	None	None		1.6 M Sodium Citrate	
<b>D 6</b>	20% w/v PEG 3350	None	None		200 mM Potassium Citrate	8.3