

## JBScreen Classic 10

Cat.-No: CS-110L

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
<b>A 1</b>	500 mM Sodium Acetate	None	100 mM Imidazole-HCl	8.0	None
<b>A 2</b>	700 mM Sodium Citrate	None	100 mM HEPES Sodium Salt	7.5	None
<b>A 3</b>	700 mM Lithium Sulfate	None	100 mM Tris-HCl	8.5	None
<b>A 4</b>	800 mM Potassium / Sodium Tartrate	None	100 mM HEPES Sodium Salt	7.5	None
<b>A 5</b>	1.0 M Ammonium dihydrogen Phosphate	None	100 mM Sodium Citrate	5.6	None
<b>A 6</b>	1.0 M Ammonium dihydrogen Phosphate	None	100 mM Tris-HCl	8.5	None
<b>B 1</b>	1.0 M Lithium Sulfate	None	100 mM Tris-HCl	8.5	10 mM Nickel (II) Chloride
<b>B 2</b>	1.0 M Sodium Acetate	None	100 mM Imidazole-HCl	8.0	None
<b>B 3</b>	1.0 M Sodium Formate	None	100 mM Sodium Acetate	4.6	None
<b>B 4</b>	1.4 M Sodium Acetate	None	100 mM MES Sodium Salt	6.5	None
<b>B 5</b>	1.4 M Sodium Citrate	None	100 mM HEPES Sodium Salt	7.5	None
<b>B 6</b>	1.5 M Lithium Sulfate	None	100 mM Tris-HCl	8.5	None
<b>C 1</b>	1.5 M Sodium Citrate	None	None	6.5	None
<b>C 2</b>	1.6 M Magnesium Sulfate	None	100 mM MES Sodium Salt	6.5	None
<b>C 3</b>	1.6 M Potassium / Sodium Tartrate	None	100 mM MES Sodium Salt	6.5	None
<b>C 4</b>	2.0 M Ammonium Formate	None	100 mM MES Sodium Salt	6.5	None
<b>C 5</b>	2.0 M Ammonium dihydrogen Phosphate	None	100 mM Tris-HCl	8.5	None
<b>C 6</b>	2.0 M Sodium Formate	None	None		None
<b>D 1</b>	2.0 M Magnesium Chloride	None	100 mM Tris-HCl	8.5	None
<b>D 2</b>	2.0 M Sodium Chloride	None	100 mM MES Sodium Salt	6.5	200 mM Sodium Acetate
<b>D 3</b>	2.0 M Sodium Formate	None	100 mM Sodium Acetate	4.6	None
<b>D 4</b>	1.0 M Ammonium dihydrogen Phosphate	30 % w/v Glycerol	100 mM Tris-HCl	8.5	None
<b>D 5</b>	4.0 M Sodium Chloride	None	100 mM HEPES Sodium Salt	7.5	None
<b>D 6</b>	3.0 M Sodium Formate	None	None		None