

JBScreen Basic 4

Cat.-No: CS-124

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
A 1	800 mM Sodium dihydrogen Phosphate	800 mM Potassium dihydrogen Phosphate	100 mM HEPES Sodium Salt	7.5	None
A 2	400 mM Ammonium dihydrogen Phosphate	None	None		None
A 3	1.0 M Ammonium dihydrogen Phosphate	None	100 mM Sodium Citrate	5.6	None
A 4	2.0 M Ammonium dihydrogen Phosphate	None	100 mM Tris-HCl	8.5	None
A 5	2.0 M Ammonium Formate	None	100 mM Sodium Acetate	4.6	None
A 6	4.0 M Ammonium Formate	None	100 mM HEPES Sodium Salt	7.5	None
B 1	2.0 M Ammonium Formate	None	None		None
B 2	500 mM Ammonium Sulfate	1.0 M Lithium Sulfate	100 mM Sodium Citrate	5.6	None
B 3	1.6 M Ammonium Sulfate	None	100 mM HEPES Sodium Salt	7.5	100 mM Sodium Chloride
B 4	1.8 M Ammonium Sulfate	None	100 mM MES Sodium Salt	6.5	10 mM Cobalt(II) Chloride
B 5	2.0 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	None
B 6	2.0 M Ammonium Sulfate	None	None		None
C 1	2.0 M Ammonium Sulfate	None	100 mM Sodium Acetate	4.6	None
C 2	2.0 M Ammonium Sulfate	None	100 mM Sodium Citrate	5.6	200 mM Sodium / Potassium Tartrate
C 3	200 mM Magnesium Formate	None	None		None
C 4	1.6 M Magnesium Sulfate	None	100 mM MES Sodium Salt	6.5	None
C 5	2.0 M Magnesium Chloride	None	100 mM Bicine	9.5	None
C 6	1.0 M Sodium Acetate	None	100 mM Imidazole-HCl	6.5	None
D 1	1.0 M Sodium Acetate	None	100 mM HEPES Sodium Salt	7.5	50 mM Cadmium Sulfate
D 2	1.4 M Sodium Acetate	None	100 mM MES Sodium Salt	6.5	None
D 3	500 mM Sodium Chloride	10 mM Magnesium Chloride	None		10 mM Hexadecyl trimethylammonium Bromide
D 4	2.0 M Sodium Chloride	None	100 mM Sodium Acetate	4.6	None
D 5	2.0 M Sodium Chloride	None	100 mM MES Sodium Salt	6.5	200 mM Sodium / Potassium dihydrogen Phosphate
D 6	4.3 M Sodium Chloride	None	100 mM HEPES Sodium Salt	7.5	None