

JBScreen Basic 1

Cat.-No: CS-121

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
A 1	25% v/v Ethylene Glycol	None	None		None
A 2	12% v/v Glycerol	1.5 M Ammonium Sulfate	100 mM Tris-HCl	8.5	None
A 3	1.0 M 1,6-Hexanediol	None	100 mM Sodium Acetate	4.6	10 mM Cobalt(II) Chloride
A 4	2.5 M 1,6-Hexanediol	None	100 mM Sodium Citrate	5.6	None
A 5	3.4 M 1,6 Hexanediol	None	100 mM Tris-HCl	8.5	200 mM Magnesium Chloride
A 6	30% v/v MPD	None	100 mM Sodium Acetate	4.6	200 mM Sodium Chloride
B 1	30% v/v MPD	None	100 mM Sodium Citrate	5.6	200 mM Ammonium Acetate
B 2	30% v/v MPD	None	100 mM Sodium Acetate	4.6	20 mM Calcium Chloride
B 3	30% v/v MPD	500 mM Ammonium Sulfate	100 mM HEPES Sodium Salt	7.5	None
B 4	30% v/v MPD	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Citrate
B 5	50% v/v MPD	None	100 mM Tris-HCl	8.5	200 mM Ammonium dihydrogen Phosphate
B 6	70 % v/v MPD	None	100 mM HEPES Sodium Salt	7.5	None
C 1	2% v/v Ethylene Imine Polymer	None	100 mM Sodium Citrate	5.6	500 mM Sodium Chloride
C 2	2% v/v PEG 400	2.0 M Ammonium Sulfate	100 mM HEPES Sodium Salt	7.5	None
C 3	28% v/v PEG 400	None	100 mM HEPES Sodium Salt	7.5	200 mM Calcium Chloride
C 4	30 % v/ v PEG 400	None	100 mM Tris-HCl	8.5	200 mM Sodium Citrate
C 5	30 % v/ v PEG 400	None	100 mM HEPES Sodium Salt	7.5	200 mM Magnesium Chloride
C 6	30 % v/ v PEG 400	None	100 mM Sodium Acetate	4.6	100 mM Calcium Chloride
D 1	20 % v/v PEG 550 MME	None	100 mM Bicine	9.5	100 mM Sodium Chloride
D 2	25% v/v PEG 550 MME	None	100 mM MES Sodium Salt	6.5	10 mM Zinc Sulfate
D 3	10% w/v PEG 1000	10% w/v PEG 8000	None		None
D 4	30% w/v PEG 1500	None	None		None
D 5	20% w/v PEG 2000 MME	None	100 mM Tris-HCl	8.5	10 mM Nickel(II) Chloride
D 6	30% w/v PEG 2000 MME	None	100 mM Sodium Acetate	4.6	200 mM Ammonium Sulfate