

JBScreen Classic 7

Cat.-No: CS-107L

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
A 1	10 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	100 mM Sodium Citrate
A 2	12 % w/v MPD	None	100 mM Tris-HCl	8.5	50 mM Magnesium Chloride
A 3	15 % w/v MPD	None	100 mM Sodium Acetate	4.6	20 mM Calcium Chloride
A 4	15 % w/v MPD	5 % w/v PEG 4000	100 mM Imidazole-HCl	8.0	None
A 5	15 % w/v MPD	None	100 mM Sodium Citrate	5.6	200 mM Ammonium Acetate
A 6	15 % w/v MPD	None	100 mM MES Sodium Salt	6.5	200 mM Magnesium Acetate
B 1	15 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Citrate
B 2	20 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	100 mM Sodium Citrate
B 3	20 % w/v MPD	None	100 mM Imidazole-HCl	8.0	None
B 4	20 % w/v MPD	4 % w/v Glycerol	None		200 mM Sodium Chloride
B 5	30 % w/v MPD	None	100 mM Sodium Acetate	4.6	20 mM Calcium Chloride
B 6	30 % w/v MPD	None	100 mM Sodium Citrate	5.6	200 mM Ammonium Acetate
C 1	30 % w/v MPD	None	100 mM MES Sodium Salt	6.5	200 mM Magnesium Acetate
C 2	30 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	500 mM Ammonium Sulfate
C 3	30 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Citrate
C 4	30 % w/v MPD	5 % w/v PEG 4000	100 mM HEPES Sodium Salt	7.5	None
C 5	30 % w/v MPD	10 % w/v PEG 4000	100 mM Imidazole-HCl	8.0	None
C 6	30 % w/v MPD	20 % w/v Ethanol	None		None
D 1	35 % w/v MPD	None	None		None
D 2	35 % w/v MPD	None	100 mM Imidazole-HCl	8.0	None
D 3	40 % w/v MPD	None	100 mM Tris-HCl	8.5	None
D 4	47 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	None
D 5	47 % w/v MPD	2 % w/v tert-Butanol	None		None
D 6	50 % w/v MPD	None	None		None