

JBScreen Classic 4

Cat.-No: CS-104L

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
A 1	25 % w/v PEG 5000 MME	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate	None
A 2	30 % w/v PEG 5000 MME	None	100 mM MES Sodium Salt	6.5	200 mM Ammonium Sulfate	None
A 3	3 % w/v PEG 6000	None	100 mM Tris-HCl	8.5	100 mM Potassium Chloride	None
A 4	10 % w/v PEG 6000	None	None		10 mM Magnesium Chloride	None
A 5	12 % w/v PEG 6000	2.0 M Sodium Chloride	None		None	None
A 6	15 % w/v PEG 6000	5 % w/v Glycerol	None		None	None
B 1	15 % w/v PEG 6000	50 mM Potassium Chloride	None		10 mM Magnesium Chloride	None
B 2	16 % w/v PEG 6000	None	None		10 mM Sodium Citrate	None
B 3	20 % w/v PEG 6000	None	50 mM Imidazole-HCl	8.0	None	None
B 4	25 % w/v PEG 6000	None	100 mM HEPES Sodium Salt	7.5	100 mM Lithium Chloride	None
B 5	28 % w/v PEG 6000	500 mM Lithium Chloride	100 mM Tris-HCl	8.5	None	None
B 6	30 % w/v PEG 6000	1.0 M Lithium Chloride	None		100 mM Sodium Acetate	None
C 1	33 % w/v PEG 6000	None	None		10 mM Sodium Citrate	None
C 2	2 % w/v PEG 8000	500 mM Lithium Sulfate	None		None	None
C 3	2 % w/v PEG 8000	1.0 M Lithium Sulfate	None		None	None
C 4	4 % w/v PEG 8000	None	None		None	None
C 5	8 % w/v PEG 8000	200 mM Lithium Chloride	None		50 mM Magnesium Sulfate	None
C 6	8 % w/v PEG 8000	None	100 mM Tris-HCl	8.5	None	None
D 1	10 % w/v PEG 8000	None	100 mM MES Sodium Salt	6.5	200 mM Zinc Acetate	None
D 2	10 % w/v PEG 8000	None	100 mM HEPES Sodium Salt	7.5	200 mM Calcium Acetate	None
D 3	10 % w/v PEG 8000	None	None		50 mM Magnesium Acetate	100 mM Sodium Acetate
D 4	10 % w/v PEG 8000	None	None		200 mM Magnesium Acetate	None
D 5	10 % w/v PEG 8000	10 % w/v Ethylene Glycol	100 mM HEPES Sodium Salt	7.5	None	None
D 6	10 % w/v PEG 8000	10 % w/v PEG 1000	None		None	None