

JBScreen Membrane 3

Cat.-No: CS-303L

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
A 1	700 mM Ammonium Sulfate	None	1.0 M Sodium / Potassium Phosphate	7.5	None	None
A 2	1.0 M Ammonium Sulfate	None	50 mM MES Sodium Salt	6.5	100 mM Zinc Acetate	None
A 3	1.2 M Ammonium Sulfate	None	50 mM Tris-HCl	7.5	None	None
A 4	1.2 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	None	None
A 5	1.4 M Ammonium Sulfate	4% w/v 2-Propanol	None		100 mM Ammonium Acetate	None
A 6	2.0 M Ammonium Sulfate	None	None		None	None
B 1	2.0 M Ammonium Sulfate	None	100 mM Sodium Citrate	5.6	None	None
B 2	2.5 M Ammonium Sulfate	2% PEG 5000 MME	100 mM HEPES Sodium Salt	7.5	None	None
B 3	3.0 M Ammonium Sulfate	None	None		None	None
B 4	3.5 M Ammonium Sulfate	None	None		None	None
B 5	3.5 M Ammonium Sulfate	None	50 mM Sodium/ Potassium Phosphate	7.5	250 mM Sodium Chloride	None
B 6	25% w/v MPD	None	100 mM Bis Tris	7.0	None	None
C 1	30% w/v MPD	None	300 mM Sodium Citrate	5.6	None	None
C 2	25% w/v Triethylene Glycol	None	None		100 mM Ammonium Sulfate	100 mM Glycine
C 3	30% w/v 2-Propanol	20% w/v Glycerol	100 mM Sodium Acetate	4.6	200 mM Calcium Chloride	None
C 4	None	None	50 mM Potassium Phosphate	8.0	None	None
C 5	None	None	100 mM Sodium Citrate	4.8	None	None
C 6	1.0 M Potassium Phosphate	None	None	6.5	1% w/v 1,4-Dioxane	None
D 1	1.0 M Sodium Citrate	None	None		None	None
D 2	1.0 M Sodium Citrate	None	None	5.6	500 mM Lithium Chloride	None
D 3	1.5 M Sodium Chloride	None	100 mM Sodium Acetate	4.6	None	None
D 4	1.5 M Potassium Phosphate	None	None	7.0	None	None
D 5	1.5 M Lithium Sulfate	None	100 mM HEPES Sodium Salt	7.5	None	None
D 6	2.0 M Sodium Chloride	None	None		100 mM Sodium Formate	None