

JBScreen Membrane 1

Cat.-No: CS-301L

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
A 1	15% w/v PEG 400	15% w/v Glycerol	100 mM HEPES Sodium Salt	7.5	200 mM Calcium Chloride	None
A 2	20% w/v PEG 400	100 mM Sodium Chloride	100 mM Sodium Citrate	5.6	20 mM Magnesium Chloride	None
A 3	25% w/v PEG 400	None	50 mM Sodium Acetate	4.6	50 mM Magnesium Acetate	None
A 4	30% w/v PEG 400	50 mM Sodium Sulfate	50 mM Tris-HCl	8.5	50 mM Lithium Sulfate	None
A 5	48% w/v PEG 400	None	100 mM HEPES Sodium Salt	7.5	200 mM Calcium Chloride	None
A 6	20% w/v PEG 550 MME	None	10 mM Tris-HCl	7.5	None	None
B 1	30% w/v PEG 550 MME	None	50 mM Tris-HCl	8.5	100 mM Magnesium Chloride	None
B 2	35% w/v PEG 600	None	None		None	None
B 3	28% w/v PEG 1000	10% w/v Glycerol	100 mM Tricine	8.0	350 mM Sodium Chloride	None
B 4	10% w/v PEG 1500	5% w/v Ethanol	None		100 mM Magnesium Chloride	100 mM Sodium Chloride
B 5	30% w/v PEG 1500	None	None		None	None
B 6	5% w/v PEG 2000	None	None		None	None
C 1	10% w/v PEG 2000	None	50 mM Tris-HCl	8.5	500 mM Magnesium Chloride	None
C 2	15% w/v PEG 2000	None	None		None	None
C 3	15% w/v PEG 2000	None	None		100 mM Lithium Chloride	None
C 4	15% w/v PEG 2000	None	100 mM Sodium Phosphate	6.2	20 mM Sodium Citrate	None
C 5	15% w/v PEG 2000	None	100 mM Sodium Phosphate	6.8	500 mM Sodium Chloride	None
C 6	15% w/v PEG 2000	None	20 mM Bis Tris	7.0	None	None
D 1	15% w/v PEG 2000	None	50 mM HEPES Sodium Salt	7.5	100 mM Magnesium Chloride	None
D 2	20% w/v PEG 2000	2% w/v MPD	100 mM Tris-HCl	8.0	300 mM Magnesium Nitrate	None
D 3	25% w/v PEG 2000	15% w/v Glycerol	100 mM Bicine	9.0	300 mM Magnesium Chloride	None
D 4	30% w/v PEG 2000	None	200 mM Sodium Phosphate	6.2	500 mM Sodium Chloride	None
D 5	8% w/v PEG 2000 MME	None	100 mM Sodium Acetate	4.6	None	None
D 6	10% w/v PEG 2000 MME	20% w/v Glycerol	100 mM Sodium Citrate	5.6	3% w/v PEG 200	None