

JBScreen Membrane HTS

Cat.-No.: CS-305L (1.7 ml per well)

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
A 7	30% w/v PEG 550 MME	None	50 mM Tris-HCl	8.5	100 mM Magnesium Chloride	None
A 8	35% w/v PEG 600	None	None		None	None
A 9	28% w/v PEG 1000	10% w/v Glycerol	100 mM Tricine	8.0	350 mM Sodium Chloride	None
A 10	10% w/v PEG 1500	5% w/v Ethanol	None		100 mM Magnesium Chloride	100 mM Sodium Chloride
A 11	30% w/v PEG 1500	None	None		None	None
A 12	5% w/v PEG 2000	None	None		None	None
B 1	10% w/v PEG 2000	None	50 mM Tris-HCl	8.5	500 mM Magnesium Chloride	None
B 2	15% w/v PEG 2000	None	None		None	None
B 3	15% w/v PEG 2000	None	None		100 mM Lithium Chloride	None
B 4	15% w/v PEG 2000	None	100 mM Sodium Phosphate	6.2	20 mM Sodium Citrate	None
B 5	15% w/v PEG 2000	None	100 mM Sodium Phosphate	6.8	500 mM Sodium Chloride	None
B 6	15% w/v PEG 2000	None	20 mM Bis Tris	7.0	None	None
B 7	15% w/v PEG 2000	None	50 mM HEPES Sodium Salt	7.5	100 mM Magnesium Chloride	None
B 8	20% w/v PEG 2000	2% w/v MPD	100 mM Tris-HCl	8.0	300 mM Magnesium Nitrate	None
B 9	25% w/v PEG 2000	15% w/v Glycerol	100 mM Bicine	9.0	300 mM Magnesium Chloride	None
B 10	30% w/v PEG 2000	None	200 mM Sodium Phosphate	6.2	500 mM Sodium Chloride	None
B 11	8% w/v PEG 2000 MME	None	100 mM Sodium Acetate	4.6	None	None
B 12	10% w/v PEG 2000 MME	20% w/v Glycerol	100 mM Sodium Citrate	5.6	3% w/v PEG 200	None

JBScreen Membrane HTS

Cat.-No.: CS-305L (1.7 ml per well)

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
C 1	12% w/v PEG 2000 MME	None	50 mM Tris-HCl	7.5	500 mM Sodium Chloride
C 2	10% w/v PEG 3350	None	50 mM Sodium Citrate	5.6	150 mM Sodium Chloride
C 3	2% w/v PEG 4000	None	50 mM Tris-HCl	7.5	None
C 4	5% w/v PEG 4000	None	None		None
C 5	5% w/v PEG 4000	None	None		100 mM Potassium Chloride
C 6	5% w/v PEG 4000	10% w/v Glycerol	50 mM MES Sodium Salt	6.5	100 mM Sodium Chloride
C 7	5% w/v PEG 4000	None	50 mM Sodium Phosphate	6.7	None
C 8	10% w/v PEG 4000	None	50 mM Tris-HCl	8.5	500 mM Sodium Chloride
C 9	12% w/v PEG 4000	None	100 mM ADA	6.5	100 mM Lithium Sulfate
C 10	12% w/v PEG 4000	None	50 mM Sodium Phosphate	6.8	None
C 11	12% w/v PEG 4000	20% w/v Glycerol	50 mM MOPS	7.0	500 mM Potassium Chloride
C 12	15% w/v PEG 4000	None	10 mM Tris-HCl	7.5	100 mM Lithium Chloride
D 1	20% w/v PEG 4000	None	100 mM Bis Tris	7.0	500 mM Sodium Chloride
D 2	20% w/v PEG 4000	None	100 mM Sodium Phosphate	7.0	500 mM Sodium Chloride
D 3	20% w/v PEG 4000	150 mM Zinc Acetate	50 mM Tris-HCl	7.5	50 mM Zinc Chloride
D 4	22% w/v PEG 4000	None	50 mM Tricine	8.0	None
D 5	22% w/v PEG 4000	None	50 mM Tris-HCl	8.5	500 mM Sodium Chloride
D 6	30% w/v PEG 4000	None	None		None
D 7	10% w/v PEG 5000 MME	None	100 mM Sodium Citrate	5.6	100 mM Magnesium Acetate
D 8	5% w/v PEG 6000	None	None		100 mM Magnesium Sulfate
D 9	10% w/v PEG 6000	150 mM Zinc Acetate	50 mM Tris-HCl	7.5	50 mM Zinc Chloride
D 10	15% w/v PEG 6000	None	50 mM Sodium Succinate	6.5	None
D 11	12% w/v PEG 8000	10% w/v MPD	None		25 mM Potassium dihydrogen Phosphate
D 12	8% w/v PEG 10000	None	100 mM Sodium Citrate	5.6	100 mM Magnesium Acetate

JBScreen Membrane HTS

Cat.-No.: CS-305L (1.7 ml per well)

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
E 1	700 mM Ammonium Sulfate	None	1.0 M Sodium / Potassium Phosphate	7.5	None	None
E 2	1.0 M Ammonium Sulfate	None	50 mM MES Sodium Salt	6.5	100 mM Zinc Acetate	None
E 3	1.2 M Ammonium Sulfate	None	50 mM Tris-HCl	7.5	None	None
E 4	1.2 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	None	None
E 5	1.4 M Ammonium Sulfate	4% w/v 2-Propanol	None		100 mM Ammonium Acetate	None
E 6	2.0 M Ammonium Sulfate	None	None		None	None
E 7	2.0 M Ammonium Sulfate	None	100 mM Sodium Citrate	5.6	None	None
E 8	2.5 M Ammonium Sulfate	2% PEG 5000 MME	100 mM HEPES Sodium Salt	7.5	None	None
E 9	3.0 M Ammonium Sulfate	None	None		None	None
E 10	3.5 M Ammonium Sulfate	None	None		None	None
E 11	3.5 M Ammonium Sulfate	None	50 mM Sodium/ Potassium Phosphate	7.5	250 mM Sodium Chloride	None
E 12	25% w/v MPD	None	100 mM Bis Tris	7.0	None	None
F 1	30% w/v MPD	None	300 mM Sodium Citrate	5.6	None	None
F 2	25% w/v Triethylene Glycol	None	None		100 mM Ammonium Sulfate	100 mM Glycine
F 3	30% w/v 2-Propanol	20% w/v Glycerol	100 mM Sodium Acetate	4.6	200 mM Calcium Chloride	None
F 4	None	None	50 mM Potassium Phosphate	8.0	None	None
F 5	None	None	100 mM Sodium Citrate	4.8	None	None
F 6	1.0 M Potassium Phosphate	None	None	6.5	1% w/v 1,4- Dioxane	None
F 7	1.0 M Sodium Citrate	None	None		None	None
F 8	1.0 M Sodium Citrate	None	None	5.6	500 mM Lithium Chloride	None
F 9	1.5 M Sodium Chloride	None	100 mM Sodium Acetate	4.6	None	None
F 10	1.5 M Potassium Phosphate	None	None	7.0	None	None
F 11	1.5 M Lithium Sulfate	None	100 mM HEPES Sodium Salt	7.5	None	None
F 12	2.0 M Sodium Chloride	None	None		100 mM Sodium Formate	None