

JBScreen JCSG++ HTS

Cat.-No.: CS-206S (1.0 ml per well)
CS-206L (1.7 ml per well)

Number	Precipitant 1	Precipitant 2	Buffer	pH	Salt	pH
A 1	50% v/v PEG 400	None	100 mM Sodium Acetate	4.5	200 mM Lithium Sulfate	
A 2	20% w/v PEG 3000	None	100 mM Sodium Citrate	5.5	None	
A 3	20% w/v PEG 3350	None	None		200 mM di-Ammonium hydrogen Citrate	5.0
A 4	30% v/v MPD	None	100 mM Sodium Acetate	4.6	20 mM Calcium Chloride	
A 5	20% w/v PEG 3350	None	None		200 mM Magnesium Formate	5.9
A 6	20% w/v PEG 1000	None	100 mM Phosphate / Citrate	4.2	200 mM Lithium Sulfate	
A 7	20% w/v PEG 8000	None	100 mM CHES	9.5	None	
A 8	20% w/v PEG 3350	None	None		200 mM Ammonium Formate	6.6
A 9	20% w/v PEG 3350	None	None		200 mM Ammonium Chloride	6.3
A 10	20% w/v PEG 3350	None	None		200 mM Potassium Formate	7.3
A 11	50% v/v MPD	None	100 mM Tris-HCl	8.5	200 mM Ammonium dihydrogen Phosphate	
A 12	20% w/v PEG 3350	None	None		200 mM Potassium Nitrate	6.9
B 1	None	None	100 mM Citric Acid	4.0	800 mM Ammonium Sulfate	
B 2	20% w/v PEG 3550	None	None		200 mM Sodium Thiocyanate	6.9
B 3	20% w/v PEG 6000	None	100 mM Bicine	9.0	None	
B 4	10% w/v PEG 8000	8% v/v Ethylene Glycol	100 mM HEPES Sodium Salt	7.5	None	
B 5	40% v/v MPD	5% w/v PEG 8000	100 mM MES Sodium Salt	6.5	None	
B 6	40% v/v Ethanol	5% w/v PEG 1000	100 mM Phosphate / Citrate	4.2	None	
B 7	8% w/v PEG 4000	None	100 mM Sodium Acetate	4.6	None	
B 8	10% w/v PEG 8000	None	100 mM Tris-HCl	7.0	200 mM Magnesium Chloride	
B 9	20% w/v PEG 6000	None	100 mM Citric Acid	5.0	None	
B 10	50% v/v PEG 200	None	100 mM MES Sodium Salt	6.5	200 mM Magnesium Chloride	
B 11	None	None	None		1.6 M Sodium Citrate	
B 12	20% w/v PEG 3350	None	None		200 mM Potassium Citrate	8.3

Data Sheet

JBScreen JCSG++ HTS

Cat.-No.: CS-206S (1.0 ml per well)

CS-206L (1.7 ml per well)

Number	Precipitant 1	Precipitant 2	Buffer	pH	Salt	pH
C 1	20% w/v PEG 8000	None	100 mM Phosphate / Citrate	4.2	200 mM Sodium Chloride	
C 2	20% w/v PEG 6000	None	100 mM Citric Acid	4.0	1.0 M Lithium Chloride	
C 3	20% w/v PEG 3350	None	None		200 mM Ammonium Nitrate	6.3
C 4	10% w/v PEG 6000	None	100 mM HEPES Sodium Salt	7.0	None	
C 5	800 mM Sodium dihydrogen Phosphate	800 mM Potassium dihydrogen Phosphate	100 mM HEPES Sodium Salt	7.5	None	
C 6	40% v/v PEG 300	None	100 mM Phosphate / Citrate	4.2	None	
C 7	10% w/v PEG 3000	None	100 mM Sodium Acetate	4.5	200 mM Zinc Acetate	
C 8	20% v/v Ethanol	None	100 mM Tris-HCl	8.5	None	
C 9	25% v/v 1,2-Propanediol	10% v/v Glycerol	100 mM Sodium / Potassium Phosphate	6.2	None	
C 10	10% w/v PEG 20000	2% v/v 1,4-Dioxane	100 mM Bicine	9.0	None	
C 11	None	None	100 mM Sodium Acetate	4.6	2.0 M Ammonium Sulfate	
C 12	10% w/v PEG 1000	10% w/v PEG 8000	None		None	
D 1	24% w/v PEG 1500	20% v/v Glycerol	None		None	
D 2	30% v/v PEG 400	None	100 mM HEPES Sodium Salt	7.5	200 mM Magnesium Chloride	
D 3	50% v/v PEG 200	None	100 mM Sodium / Potassium Phosphate	6.2	200 mM Sodium Chloride	
D 4	30% w/v PEG 8000	None	100 mM Sodium Acetate	4.5	200 mM Lithium Sulfate	
D 5	70% v/v MPD	None	100 mM HEPES Sodium Salt	7.5	None	
D 6	20% w/v PEG 8000	None	100 mM Tris-HCl	8.5	200 mM Magnesium Chloride	
D 7	40% v/v PEG 400	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate	
D 8	40% v/v MPD	None	100 mM Tris-HCl	8.0	None	
D 9	25.5% w/v PEG 4000	15% v/v Glycerol	None		170 mM Ammonium Sulfate	
D 10	40% v/v PEG 300	None	100 mM MES Sodium Salt	6.5	200 mM Calcium Acetate	
D 11	14% v/v 2-Propanol	30% v/v Glycerol	70 mM Sodium Acetate	4.6	140 mM Calcium Chloride	
D 12	16% w/v PEG 8000	20% v/v Glycerol	None		40 mM Potassium dihydrogen Phosphate	

JBScreen JCSG++ HTS

Cat.-No.: CS-206S (1.0 ml per well)

CS-206L (1.7 ml per well)

Number	Precipitant 1	Precipitant 2	Buffer	pH	Salt	pH
E 1	None	None	100 mM MES Sodium Salt	6.5	1.0 M Sodium Citrate	
E 2	2.0 M Ammonium Sulfate	None	100 mM MES Sodium Salt	6.5	200 mM Sodium Chloride	
E 3	10% v/v 2-Propanol	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Chloride	
E 4	1.26 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	200 mM Lithium Sulfate	
E 5	40% v/v MPD	None	100 mM CAPS	10.5	None	
E 6	20% w/v PEG 3000	None	100 mM Imidazole-HCl	8.0	200 mM Zinc Acetate	
E 7	10% v/v 2-Propanol	None	100 mM MES Sodium Salt	6.5	200 mM Zinc Acetate	
E 8	None	None	100 mM Sodium Acetate	4.5	1.0 M di-Ammonium hydrogen Phosphate	
E 9	None	None	100 mM MES Sodium Salt	6.5	1.6 M Magnesium Sulfate	
E 10	10% w/v PEG 6000	None	100 mM Bicine	9.0	None	
E 11	14.4% w/v PEG 8000	20% v/v Glycerol	80 mM MES Sodium Salt	6.5	160 mM Calcium Acetate	
E 12	10% w/v PEG 8000	None	100 mM Imidazole-HCl	8.0	None	
F 1	30% v/v Jeffamine M-600	None	100 mM MES Sodium Salt	6.5	50 mM Cesium Chloride	
F 2	None	None	100 mM Citric Acid	5.0	3.15 M Ammonium Sulfate	
F 3	20% v/v MPD	None	100 mM Tris-HCl	8.0	None	
F 4	20% v/v Jeffamine M-600	None	100 mM HEPES Sodium Salt	7.5	None	
F 5	50% v/v Ethylene Glycol	None	100 mM Tris-HCl	8.5	200 mM Magnesium Chloride	
F 6	10% v/v MPD	None	100 mM Bicine	9.0	None	
F 7	None	None	None		800 mM Succinic Acid	7.0
F 8	None	None	None		2.1 M D,L-Malic Acid	7.0
F 9	None	None	None		2.4 M Sodium Malonate	7.0
F 10	0.5% w/v Jeffamine ED-2001 pH 7.0	None	100 mM HEPES Sodium Salt	7.0	1.1 M Sodium Malonate	7.0
F 11	1% w/v PEG 2000 MME	None	100 mM HEPES Sodium Salt	7.0	1.0 M Succinic Acid	7.0
F 12	30% v/v Jeffamine M-600 pH 7.0	None	100 mM HEPES Sodium Salt	7.0	None	

JBScreen JCSG++ HTS

Cat.-No.: CS-206S (1.0 ml per well)

CS-206L (1.7 ml per well)

Number	Precipitant 1	Buffer	pH	Salt	pH
G 1	30% w/v Jeffamine ED-2001 pH 7.0	100 mM HEPES Sodium Salt	7.0	None	
G 2	22% w/v Polyacrylic Acid 5100 Sodium Salt	100 mM HEPES Sodium Salt	7.5	200 mM Magnesium Chloride	
G 3	20% w/v Polyvinylpyrrolidone K 15	100 mM Tris-HCl	8.5	100 mM Cobalt Chloride	
G 4	20% w/v PEG 2000 MME	100 mM Tris-HCl	8.5	200 mM Trimethyl-N-oxide	
G 5	12% w/v PEG 3350	100 mM HEPES Sodium Salt	7.5	5 mM Cobalt Chloride, 5 mM Cadmium Chloride 5 mM Magnesium Chloride, 5 mM Nickel Chloride	
G 6	20% w/v PEG 3350	None		200 mM Sodium Malonate	7.0
G 7	15% w/v PEG 3350	None		100 mM Succinic Acid	7.0
G 8	20% w/v PEG 3350	None		150 mM D,L-Malic Acid	7.0
G 9	30% w/v PEG 2000 MME	None		100 mM Potassium Thiocyanate	
G 10	30% w/v PEG 2000 MME	None		150 mM Potassium Bromide	
G 11	None	100 mM Bis-Tris	5.5	2.0 M Ammonium Sulfate	
G 12	None	100 mM Bis-Tris	5.5	3.0 M Sodium Chloride	
H 1	None	100 mM Bis-Tris	5.5	300 mM Magnesium Formate	
H 2	1% w/v PEG 3550	100 mM Bis-Tris	5.5	1.0 M Ammonium Sulfate	
H 3	25% w/v PEG 3350	100 mM Bis-Tris	5.5	None	
H 4	45% v/v MPD	100 mM Bis-Tris	5.5	200 mM Calcium Chloride	
H 5	45% v/v MPD	100 mM Bis-Tris	5.5	200 mM Ammonium Acetate	
H 6	17% w/v PEG 10000	100 mM Bis-Tris	5.5	100 mM Ammonium Acetate	
H 7	25% w/v PEG 3350	100 mM Bis-Tris	5.5	200 mM Ammonium Sulfate	
H 8	25% w/v PEG 3350	100 mM Bis-Tris	5.5	200 mM Sodium Chloride	
H 9	25% w/v PEG 3350	100 mM Bis-Tris	5.5	200 mM Lithium Sulfate	
H 10	25% w/v PEG 3350	100 mM Bis-Tris	5.5	200 mM Ammonium Acetate	
H 11	25% w/v PEG 3350	100 mM Bis-Tris	5.5	200 mM Magnesium Chloride	
H 12	45% v/v MPD	100 mM HEPES	7.5	200 mM Ammonium Acetate	