

JBScreen PACT++ HTS

Cat.-No.: CS-207S (1.0 ml per well)
CS-207L (1.7 ml per well)

Number	Precipitant	Buffer	pH	Salt
A 1	25% w/v PEG 1500	100 mM SPG Buffer	4.0	None
A 2	25% w/v PEG 1500	100 mM SPG Buffer	5.0	None
A 3	25% w/v PEG 1500	100 mM SPG Buffer	6.0	None
A 4	25% w/v PEG 1500	100 mM SPG Buffer	7.0	None
A 5	25% w/v PEG 1500	100 mM SPG Buffer	8.0	None
A 6	25% w/v PEG 1500	100 mM SPG Buffer	9.0	None
A 7	20% w/v PEG 6000	100 mM Sodium Acetate	5.0	200 mM Sodium Chloride
A 8	20% w/v PEG 6000	100 mM Sodium Acetate	5.0	200 mM Ammonium Chloride
A 9	20% w/v PEG 6000	100 mM Sodium Acetate	5.0	200 mM Lithium Chloride
A 10	20% w/v PEG 6000	100 mM Sodium Acetate	5.0	200 mM Magnesium Chloride
A 11	20% w/v PEG 6000	100 mM Sodium Acetate	5.0	200 mM Calcium Chloride
A 12	20% w/v PEG 6000	100 mM Sodium Acetate	5.0	10 mM Zinc Chloride
B 1	25% w/v PEG 1500	100 mM MIB Buffer	4.0	None
B 2	25% w/v PEG 1500	100 mM MIB Buffer	5.0	None
B 3	25% w/v PEG 1500	100 mM MIB Buffer	6.0	None
B 4	25% w/v PEG 1500	100 mM MIB Buffer	7.0	None
B 5	25% w/v PEG 1500	100 mM MIB Buffer	8.0	None
B 6	25% w/v PEG 1500	100 mM MIB Buffer	9.0	None
B 7	20% w/v PEG 6000	100 mM MES Sodium Salt	6.0	200 mM Sodium Chloride
B 8	20% w/v PEG 6000	100 mM MES Sodium Salt	6.0	200 mM Ammonium Chloride
B 9	20% w/v PEG 6000	100 mM MES Sodium Salt	6.0	200 mM Lithium Chloride
B 10	20% w/v PEG 6000	100 mM MES Sodium Salt	6.0	200 mM Magnesium Chloride
B 11	20% w/v PEG 6000	100 mM MES Sodium Salt	6.0	200 mM Calcium Chloride
B 12	20% w/v PEG 6000	100 mM MES Sodium Salt	6.0	10 mM Zinc Chloride

JBScreen PACT++ HTS

Cat.-No.: CS-207S (1.0 ml per well)
CS-207L (1.7 ml per well)

Number	Precipitant	Buffer	pH	Salt
C 1	25% w/v PEG 1500	100 mM TBG Buffer	4.0	None
C 2	25% w/v PEG 1500	100 mM TBG Buffer	5.0	None
C 3	25% w/v PEG 1500	100 mM TBG Buffer	6.0	None
C 4	25% w/v PEG 1500	100 mM TBG Buffer	7.0	None
C 5	25% w/v PEG 1500	100 mM TBG Buffer	8.0	None
C 6	25% w/v PEG 1500	100 mM TBG Buffer	9.0	None
C 7	20% w/v PEG 6000	100 mM HEPES Sodium Salt	7.0	200 mM Sodium Chloride
C 8	20% w/v PEG 6000	100 mM HEPES Sodium Salt	7.0	200 mM Ammonium Chloride
C 9	20% w/v PEG 6000	100 mM HEPES Sodium Salt	7.0	200 mM Lithium Chloride
C 10	20% w/v PEG 6000	100 mM HEPES Sodium Salt	7.0	200 mM Magnesium Chloride
C 11	20% w/v PEG 6000	100 mM HEPES Sodium Salt	7.0	200 mM Calcium Chloride
C 12	20% w/v PEG 6000	100 mM HEPES Sodium Salt	7.0	10 mM Zinc Chloride
D 1	25% w/v PEG 1500	100 mM MMT Buffer	4.0	None
D 2	25% w/v PEG 1500	100 mM MMT Buffer	5.0	None
D 3	25% w/v PEG 1500	100 mM MMT Buffer	6.0	None
D 4	25% w/v PEG 1500	100 mM MMT Buffer	7.0	None
D 5	25% w/v PEG 1500	100 mM MMT Buffer	8.0	None
D 6	25% w/v PEG 1500	100 mM MMT Buffer	9.0	None
D 7	20% w/v PEG 6000	100 mM Tris HCl	8.0	200 mM Sodium Chloride
D 8	20% w/v PEG 6000	100 mM Tris HCl	8.0	200 mM Ammonium Chloride
D 9	20% w/v PEG 6000	100 mM Tris HCl	8.0	200 mM Lithium Chloride
D 10	20% w/v PEG 6000	100 mM Tris HCl	8.0	200 mM Magnesium Chloride
D 11	20% w/v PEG 6000	100 mM Tris HCl	8.0	200 mM Calcium Chloride
D 12	20% w/v PEG 6000	100 mM Tris HCl	8.0	10 mM Zinc Chloride

JBScreen PACT++ HTS

Cat.-No.: CS-207S (1.0 ml per well)
CS-207L (1.7 ml per well)

Number	Precipitant	Buffer	pH	Salt
E 1	20% w/v PEG 3350	None		200 mM Sodium Fluoride
E 2	20% w/v PEG 3350	None		200 mM Sodium Bromide
E 3	20% w/v PEG 3350	None		200 mM Sodium Iodide
E 4	20% w/v PEG 3350	None		200 mM Potassium Thiocyanate
E 5	20% w/v PEG 3350	None		200 mM Sodium Nitrate
E 6	20% w/v PEG 3350	None		200 mM Sodium Formate
E 7	20% w/v PEG 3350	None		200 mM Sodium Acetate
E 8	20% w/v PEG 3350	None		200 mM Sodium Sulfate
E 9	20% w/v PEG 3350	None		200 mM Potassium/Sodium Tartrate
E 10	20% w/v PEG 3350	None		200 mM Sodium/Potassium Phosphate
E 11	20% w/v PEG 3350	None		200 mM Sodium Citrate
E 12	20% w/v PEG 3350	None		200 mM Sodium Malonate
F 1	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Fluoride
F 2	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Bromide
F 3	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Iodide
F 4	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Potassium Thiocyanate
F 5	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Nitrate
F 6	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Formate
F 7	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Acetate
F 8	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Sulfate
F 9	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Potassium/Sodium Tartrate
F 10	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium/Potassium Phosphate
F 11	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Citrate
F 12	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Malonate

JBScreen PACT++ HTS

Cat.-No.: CS-207S (1.0 ml per well)
CS-207L (1.7 ml per well)

Number	Precipitant	Buffer	pH	Salt
E 1	20% w/v PEG 3350	None		200 mM Sodium Fluoride
E 2	20% w/v PEG 3350	None		200 mM Sodium Bromide
E 3	20% w/v PEG 3350	None		200 mM Sodium Iodide
E 4	20% w/v PEG 3350	None		200 mM Potassium Thiocyanate
E 5	20% w/v PEG 3350	None		200 mM Sodium Nitrate
E 6	20% w/v PEG 3350	None		200 mM Sodium Formate
E 7	20% w/v PEG 3350	None		200 mM Sodium Acetate
E 8	20% w/v PEG 3350	None		200 mM Sodium Sulfate
E 9	20% w/v PEG 3350	None		200 mM Potassium/Sodium Tartrate
E 10	20% w/v PEG 3350	None		200 mM Sodium/Potassium Phosphate
E 11	20% w/v PEG 3350	None		200 mM Sodium Citrate
E 12	20% w/v PEG 3350	None		200 mM Sodium Malonate
F 1	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Fluoride
F 2	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Bromide
F 3	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Iodide
F 4	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Potassium Thiocyanate
F 5	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Nitrate
F 6	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Formate
F 7	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Acetate
F 8	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Sulfate
F 9	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Potassium/Sodium Tartrate
F 10	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium/Potassium Phosphate
F 11	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Citrate
F 12	20% w/v PEG 3350	100 mM Bis-Tris Propane	6.5	200 mM Sodium Malonate