

JBScreen PACT++ 2

Cat.-No: CS-162

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