

JBScreen Basic 3

Cat.-No: CS-123

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
A 1	10% w/v PEG 10000	2 %v/v 1,4-Dioxane	100 mM Bicine	9.5	None
A 2	20% w/v PEG 10000	None	100 mM HEPES Sodium Salt	7.5	None
A 3	12% w/v PEG 20000	None	100 mM MES Sodium Salt	6.5	None
A 4	5% v/v 2-Propanol	2.0 M Ammonium Sulfate	None		None
A 5	20% v/v 2-Propanol	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Citrate
A 6	20% v/v 2-Propanol	None	100 mM Sodium Acetate	4.6	200 mM Calcium Chloride
B 1	30% v/v 2-Propanol	None	100 mM HEPES Sodium Salt	7.5	200 mM Magnesium Chloride
B 2	30% v/v 2-Propanol	None	100 mM Tris-HCl	8.5	200 mM Ammonium Acetate
B 3	10% v/v 1,4-Dioxane	1.6 M Ammonium Sulfate	100 mM MES Sodium Salt	6.5	None
B 4	35% v/v 1,4-Dioxane	None	None		None
B 5	10% v/v Ethanol	1.5 M Sodium Chloride	None		None
B 6	20% v/v Ethanol	None	100 mM Tris-HCl	8.5	None
C 1	25% v/v tert-Butanol	None	100 mM Tris-HCl	8.5	None
C 2	35% v/v tert-Butanol	None	100 mM Sodium Citrate	5.6	None
C 3	1.0 M Imidazole-HCl	None	None	7.0	None
C 4	1.0 M Lithium Sulfate	None	100 mM Tris-HCl	8.5	10 mM Nickel(II)-Chloride
C 5	1.5 M Lithium Sulfate	None	100 mM HEPES Sodium Salt	7.5	None
C 6	400 mM Sodium / Potassium Tartrate	None	None		None
D 1	800mM Sodium / Potassium Tartrate	None	100 mM HEPES Sodium Salt	7.5	None
D 2	1.4 M Sodium Citrate	None	100 mM HEPES Sodium Salt	7.5	None
D 3	1.6 M Sodium Citrate	None	None	6.5	None
D 4	10% v/v Jeffamine M-600	None	100 mM Sodium Citrate	5.6	10 mM Ferric(III)-Chloride
D 5	20% v/v Jeffamine M-600	None	100 mM HEPES Sodium Salt	7.5	None
D 6	30% v/v Jeffamine M-600	None	100 mM MES Sodium Salt	6.5	50 mM Cesium Chloride