



## JBScreen Basic 4

Cat.-No.: CS-124

# SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
G1	800 mM Potassium di-hydrogen phosphate	800 mM Sodium di-hydrogen phosphate	100 mM HEPES; pH 7.5	none
G2	400 mM Ammonium di-hydrogen phosphate	none	none	none
G3	1 M Ammonium di-hydrogen phosphate	none	100 mM tri-Sodium citrate; pH 5.6	none
G4	2 M Ammonium di-hydrogen phosphate	none	100 mM TRIS; pH 8.5	none
G5	2 M Ammonium formate	none	100 mM Sodium acetate; pH 4.6	none
G6	4 M Ammonium formate	none	100 mM HEPES; pH 7.5	none
G7	2 M Ammonium formate	none	none	none
G8	500 mM Ammonium sulfate	1 M Lithium sulfate	100 mM tri-Sodium citrate; pH 5.6	none
G9	1.6 M Ammonium sulfate	none	100 mM HEPES; pH 7.5	100 mM Sodium chloride
G10	1.8 M Ammonium sulfate	none	100 mM MES; pH 6.5	10 mM Cobalt (II) chloride
G11	2 M Ammonium sulfate	none	100 mM TRIS; pH 8.5	none
G12	2 M Ammonium sulfate	none	none	none
H1	2 M Ammonium sulfate	none	100 mM Sodium acetate; pH 4.6	none
H2	2 M Ammonium sulfate	none	100 mM tri-Sodium citrate; pH 5.6	200 mM Potassium Sodium tartrate
H3	200 mM Magnesium formate	none	none	none
H4	1.6 M Magnesium sulfate	none	100 mM MES; pH 6.5	none
H5	2 M Magnesium chloride	none	100 mM BICINE; pH 9.5	none
H6	1 M Sodium acetate	none	100 mM Imidazole; pH 6.5	none
H7	1 M Sodium acetate	none	100 mM HEPES; pH 7.5	50 mM Cadmium sulfate
H8	1.4 M Sodium acetate	none	100 mM MES; pH 6.5	none
H9	500 mM Sodium chloride	10 mM Magnesium chloride	none	10 mM Cetyltrimethylammonium bromide
H10	2 M Sodium chloride	none	100 mM Sodium acetate; pH 4.6	none
H11	2 M Sodium chloride	none	100 mM MES; pH 6.5	100 mM Sodium di-hydrogen phosphate, 100 mM Potassium di-hydrogen phosphate
H12	4.3 M Sodium chloride	none	100 mM HEPES; pH 7.5	none

\*pH values indicated are those of the 1.0 M buffer stock solution prior to dilution with other components

