



JBScreen Nuc-Pro 4

Cat.-No.: CS-184

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
G1	10 % w/v 1,6-Hexanediol	none	50 mM MES; pH 6.5	20 mM Magnesium chloride
G2	20 % w/v 1,6-Hexanediol	none	50 mM HEPES; pH 7.0	50 mM Ammonium chloride, 10 mM Magnesium chloride
G3	35 % w/v 1,6-Hexanediol	none	50 mM TRIS; pH 8.5	75 mM Magnesium sulfate
G4	10 % v/v 2-Methyl-2,4-pentanediol	none	50 mM TRIS; pH 7.5	50 mM Ammonium acetate
G5	10 % v/v 2-Methyl-2,4-pentanediol	none	50 mM HEPES; pH 7.0	80 mM Potassium chloride, 10 mM Magnesium sulfate
G6	15 % v/v 2-Methyl-2,4-pentanediol	none	50 mM ADA; pH 6.5	100 mM Sodium acetate
G7	15 % v/v 2-Methyl-2,4-pentanediol	none	50 mM di-Sodium succinate; pH 5.5	10 mM Magnesium acetate
G8	18 % v/v 2-Methyl-2,4-pentanediol	none	20 mM MES; pH 5.8	10 mM Magnesium chloride
G9	23 % v/v 2-Methyl-2,4-pentanediol	none	50 mM MES; pH 6.0	100 mM Sodium chloride
G10	26 % v/v 2-Methyl-2,4-pentanediol	none	50 mM PIPES; pH 7.0	65 mM Magnesium chloride, 1 mM Cobalt (III) Hexamine chloride
G11	27 % v/v 2-Methyl-2,4-pentanediol	none	20 mM MES; pH 5.8	400 mM Sodium chloride, 120 mM Calcium chloride
G12	35 % v/v 2-Methyl-2,4-pentanediol	none	20 mM BIS-TRIS; pH 6.0	50 mM Sodium chloride, 10 mM Calcium chloride
H1	50 % v/v 2-Methyl-2,4-pentanediol	none	100 mM tri-Sodium citrate; pH 5.6	10 mM Magnesium chloride
H2	5 % v/v 2-Propanol	none	50 mM TRIS; pH 7.5	10 mM Magnesium chloride
H3	5 % v/v 2-Propanol	none	50 mM MES; pH 6.5	100 mM Calcium acetate
H4	9 % v/v 2-Propanol	none	50 mM Imidazole; pH 7.2	15 mM Magnesium acetate, 15 mM Magnesium chloride
H5	10 % v/v 2-Propanol	none	50 mM MES; pH 6.5	80 mM Ammonium acetate
H6	10 % v/v 2-Propanol	none	50 mM di-Sodium succinate; pH 5.5	2 mM Cobalt (II) chloride
H7	13 % v/v 2-Propanol	none	50 mM MOPS; pH 7.0	200 mM Potassium chloride, 6 mM Cobalt (III) Hexamine chloride
H8	15 % v/v 2-Propanol	none	50 mM MES; pH 6.0	20 mM Magnesium chloride
H9	10 % v/v 1,4-Dioxane	none	50 mM HEPES; pH 7.5	none
H10	10 % v/v Ethanol	none	50 mM MES; pH 6.5	20 mM Magnesium chloride, 1 mM Cobalt (II) chloride
H11	20 % v/v Ethylene glycol	5 % w/v Polyethylene glycol 3,350	none	20 mM Magnesium chloride
H12	15 % v/v Glycerol	none	100 mM Sodium acetate; pH 4.6	200 mM Sodium chloride

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

