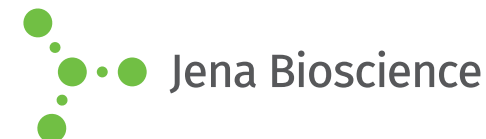




JBScreen Nuc-Pro 2

Cat.-No.: CS-182

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
C1	20 % w/v Polyethylene glycol 4,000	100 mM tri-Sodium citrate; pH 5.6	100 mM Potassium chloride, 10 % v/v 2-Propanol
C2	20 % w/v Polyethylene glycol 4,000	50 mM MOPS; pH 7.0	100 mM Sodium chloride
C3	20 % w/v Polyethylene glycol 4,000	100 mM HEPES; pH 7.5	none
C4	24 % w/v Polyethylene glycol 4,000	50 mM di-Sodium succinate; pH 5.5	60 mM Magnesium chloride, 300 mM Sodium chloride
C5	25 % w/v Polyethylene glycol 4,000	50 mM tri-Sodium citrate; pH 5.0	50 mM Ammonium sulfate
C6	30 % w/v Polyethylene glycol 4,000	50 mM MES; pH 6.5	80 mM Magnesium acetate
C7	30 % w/v Polyethylene glycol 4,000	50 mM TRIS; pH 8.5	150 mM Ammonium chloride, 10 mM Calcium chloride
C8	32 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	5 % v/v Glycerol
C9	36 % w/v Polyethylene glycol 4,000	50 mM Sodium acetate; pH 5.0	none
C10	5 % w/v Polyethylene glycol 6,000	20 mM BIS-TRIS; pH 6.0	60 mM di-Ammonium hydrogen citrate, 20 mM Magnesium chloride
C11	10 % w/v Polyethylene glycol 6,000	50 mM tri-Sodium citrate; pH 4.0	100 mM Sodium chloride
C12	10 % w/v Polyethylene glycol 6,000	50 mM HEPES; pH 7.0	200 mM Ammonium acetate, 150 mM Magnesium acetate
D1	15 % w/v Polyethylene glycol 6,000	10 mM TRIS; pH 7.5	none
D2	18 % w/v Polyethylene glycol 6,000	50 mM Sodium acetate; pH 5.0	none
D3	20 % w/v Polyethylene glycol 6,000	50 mM BIS-TRIS propane; pH 7.0	7 % v/v 2-Methyl-2,4-pentanediol, 5 % v/v 2-Methyl-2-propanol
D4	20 % w/v Polyethylene glycol 6,000	50 mM di-Potassium L-malate; pH 5.0	30 mM Calcium chloride
D5	5 % w/v Polyethylene glycol 8,000	50 mM HEPES; pH 7.5	20 mM Magnesium chloride
D6	10 % w/v Polyethylene glycol 8,000	100 mM TRIS; pH 8.0	10 % v/v Glycerol, 1 mM Tris(2-carboxyethyl)phosphine hydrochloride
D7	10 % w/v Polyethylene glycol 8,000	50 mM MES; pH 6.5	200 mM Potassium chloride, 100 mM Magnesium acetate
D8	15 % w/v Polyethylene glycol 8,000	50 mM BIS-TRIS propane; pH 6.8	100 mM Ammonium sulfate, 10 % v/v Glycerol
D9	15 % w/v Polyethylene glycol 8,000	100 mM MES; pH 6.5	200 mM Calcium acetate
D10	18 % w/v Polyethylene glycol 8,000	100 mM TRIS; pH 8.0	200 mM Magnesium formate
D11	20 % w/v Polyethylene glycol 8,000	10 mM TRIS; pH 7.5	10 mM Calcium chloride
D12	15 % w/v Polyethylene glycol 20,000	100 mM MES; pH 6.5	60 mM Manganese (II) chloride

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

