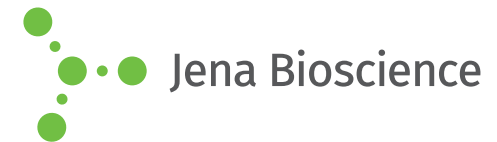




JBScreen JCSG++ 1

Cat.-No.: CS-151

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
A1	50 % v/v Polyethylene glycol 400	none	100 mM Sodium acetate; pH 4.5	200 mM Lithium sulfate
A2	20 % w/v Polyethylene glycol 3,000	none	100 mM tri-Sodium citrate; pH 5.5	none
A3	20 % w/v Polyethylene glycol 3,350	none	none	200 mM di-Ammonium hydrogen citrate; pH 5.0
A4	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM Sodium acetate; pH 4.6	20 mM Calcium chloride
A5	20 % w/v Polyethylene glycol 3,350	none	none	200 mM Magnesium formate; pH 5.9
A6	20 % w/v Polyethylene glycol 1,000	none	100 mM Potassium phosphate citrate; pH 4.2	200 mM Lithium sulfate
A7	20 % w/v Polyethylene glycol 8,000	none	100 mM CHES; pH 9.5	none
A8	20 % w/v Polyethylene glycol 3,350	none	none	200 mM Ammonium formate; pH 6.6
A9	20 % w/v Polyethylene glycol 3,350	none	none	200 mM Ammonium chloride; pH 6.3
A10	20 % w/v Polyethylene glycol 3,350	none	none	200 mM Potassium formate; pH 7.3
A11	50 % v/v 2-Methyl-2,4-pentanediol	none	100 mM TRIS; pH 8.5	200 mM Ammonium di-hydrogen phosphate
A12	20 % w/v Polyethylene glycol 3,350	none	none	200 mM Potassium nitrate; pH 6.9
B1	none	none	100 mM tri-Sodium citrate; pH 4.0	800 mM Ammonium sulfate
B2	20 % w/v Polyethylene glycol 3,350	none	none	200 mM Sodium thiocyanate; pH 6.9
B3	20 % w/v Polyethylene glycol 6,000	none	100 mM BICINE; pH 9.0	none
B4	10 % w/v Polyethylene glycol 8,000	8 % v/v Ethylene glycol	100 mM HEPES; pH 7.5	none
B5	40 % v/v 2-Methyl-2,4-pentanediol	5 % w/v Polyethylene glycol 8,000	100 mM MES; pH 6.5	none
B6	30 % v/v Ethanol	5 % w/v Polyethylene glycol 1,000	100 mM Potassium phosphate citrate; pH 4.2	none
B7	8 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	none
B8	10 % w/v Polyethylene glycol 8,000	none	100 mM TRIS; pH 7.0	200 mM Magnesium chloride
B9	20 % w/v Polyethylene glycol 6,000	none	100 mM tri-Sodium citrate; pH 5.0	none
B10	50 % v/v Polyethylene glycol 200	none	100 mM MES; pH 6.5	200 mM Magnesium chloride
B11	none	none	none	1.6 M tri-Sodium citrate



JBScreen JCSG++ 1

Cat.-No.: CS-151

SCREEN FORMULATION



B12	20 % w/v Polyethylene glycol 3,350	none	none	200 mM tri-Potassium citrate; pH 8.3
-----	------------------------------------	------	------	--------------------------------------

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