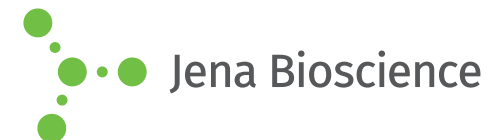




## SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
A1	25 % v/v Ethylene glycol	none	none	none
A2	12 % v/v Glycerol	1.5 M Ammonium sulfate	100 mM TRIS; pH 8.5	none
A3	1 M 1,6-Hexanediol	none	100 mM Sodium acetate; pH 4.6	10 mM Cobalt (II) chloride
A4	2.5 M 1,6-Hexanediol	none	100 mM tri-Sodium citrate; pH 5.6	none
A5	3.4 M 1,6-Hexanediol	none	100 mM TRIS; pH 8.5	200 mM Magnesium chloride
A6	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM Sodium acetate; pH 4.6	200 mM Sodium chloride
A7	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium acetate
A8	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM Sodium acetate; pH 4.6	20 mM Calcium chloride
A9	30 % v/v 2-Methyl-2,4-pentanediol	500 mM Ammonium sulfate	100 mM HEPES; pH 7.5	none
A10	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM HEPES; pH 7.5	200 mM tri-Sodium citrate
A11	50 % v/v 2-Methyl-2,4-pentanediol	none	100 mM TRIS; pH 8.5	200 mM Ammonium di-hydrogen phosphate
A12	70 % v/v 2-Methyl-2,4-pentanediol	none	100 mM HEPES; pH 7.5	none
B1	2 % w/v Ethylene imine polymer	none	100 mM tri-Sodium citrate; pH 5.6	500 mM Sodium chloride
B2	2 % v/v Polyethylene glycol 400	2 M Ammonium sulfate	100 mM HEPES; pH 7.5	none
B3	28 % v/v Polyethylene glycol 400	none	100 mM HEPES; pH 7.5	200 mM Calcium chloride
B4	30 % v/v Polyethylene glycol 400	none	100 mM TRIS; pH 8.5	200 mM tri-Sodium citrate
B5	30 % v/v Polyethylene glycol 400	none	100 mM HEPES; pH 7.5	200 mM Magnesium chloride
B6	30 % v/v Polyethylene glycol 400	none	100 mM Sodium acetate; pH 4.6	100 mM Calcium chloride
B7	20 % v/v Polyethylene glycol monomethyl ether 550	none	100 mM BICINE; pH 9.5	100 mM Sodium chloride
B8	25 % v/v Polyethylene glycol monomethyl ether 550	none	100 mM MES; pH 6.5	10 mM Zinc sulfate
B9	10 % w/v Polyethylene glycol 1,000	10 % w/v Polyethylene glycol 8,000	none	none
B10	30 % w/v Polyethylene glycol 1,500	none	none	none
B11	20 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM TRIS; pH 8.5	10 mM Nickel (II) chloride
B12	30 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate

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