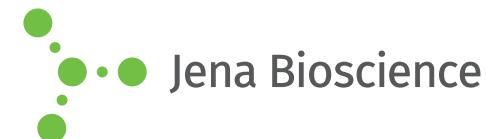




JBScreen Nuc-Pro 3

Cat.-No.: CS-183

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
E1	1.2 M Ammonium sulfate	none	100 mM HEPES; pH 7.5	2 % w/v Polyethylene glycol 400
E2	20 mM Calcium chloride	none	100 mM Glycine; pH 8.0	none
E3	2 M Lithium chloride	none	50 mM MES; pH 6.0	200 mM Calcium acetate, 1 mM Cobalt (II) chloride
E4	600 mM Lithium sulfate	none	50 mM MES; pH 6.0	10 mM Magnesium chloride
E5	1 M Lithium sulfate	8 % w/v Polyethylene glycol 400	50 mM HEPES; pH 7.5	none
E6	1 M Lithium sulfate	none	none	50 mM tri-Sodium citrate, 3 % w/v 2-Propanol
E7	1.2 M Lithium sulfate	none	50 mM MES; pH 6.5	30 mM Magnesium chloride
E8	1.2 M Lithium sulfate	none	50 mM MES; pH 6.5	50 mM Magnesium chloride, 2 mM Cobalt (II) chloride
E9	1.5 M Lithium sulfate	none	50 mM TRIS; pH 8.5	5 % w/v Glycerol
E10	1.6 M Lithium sulfate	2 % w/v Polyethylene glycol 1,000	50 mM HEPES; pH 7.5	none
E11	1.7 M Lithium sulfate	none	50 mM HEPES; pH 7.0	50 mM Magnesium sulfate
E12	1.7 M Lithium sulfate	none	none	10 % w/v Glycerol
F1	2 M Lithium sulfate	none	none	3 % w/v 2-Methyl-2,4-pentanediol
F2	5 mM Magnesium chloride	none	50 mM MES; pH 6.5	2.5 mM Cobalt (II) chloride
F3	40 mM Magnesium chloride	none	50 mM MES; pH 6.0	none
F4	80 mM Magnesium chloride	none	50 mM HEPES; pH 7.5	none
F5	200 mM Sodium chloride	none	50 mM BIS-TRIS propane; pH 7.0	none
F6	400 mM Sodium chloride	none	50 mM BIS-TRIS propane; pH 6.8	none
F7	600 mM Sodium chloride	none	50 mM MES; pH 6.0	100 mM Ammonium acetate, 5 mM Magnesium sulfate
F8	2.5 M Sodium chloride	none	50 mM TRIS; pH 7.5	200 mM Magnesium chloride
F9	1.8 M Sodium formate	none	100 mM TRIS; pH 8.0	none
F10	2 M Sodium formate	none	100 mM Sodium acetate; pH 4.6	none
F11	100 mM di-Sodium hydrogen phosphate	none	none	80 mM Sodium chloride
F12	1 M di-Sodium tartrate	none	50 mM TRIS; pH 7.5	30 mM Magnesium chloride

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

