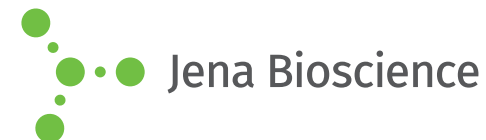




## JBScreen JCSG++ 3

Cat.-No.: CS-153

# SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
E1	none	none	100 mM MES; pH 6.5	1 M tri-Sodium citrate
E2	2 M Ammonium sulfate	none	100 mM MES; pH 6.5	200 mM Sodium chloride
E3	10 % v/v 2-Propanol	none	100 mM HEPES; pH 7.5	200 mM Sodium chloride
E4	1.26 M Ammonium sulfate	none	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
E5	40 % v/v 2-Methyl-2,4-pentanediol	none	100 mM CAPS; pH 10.5	none
E6	20 % w/v Polyethylene glycol 3,000	none	100 mM Imidazole; pH 8.0	200 mM Zinc acetate
E7	10 % v/v 2-Propanol	none	100 mM MES; pH 6.5	200 mM Zinc acetate
E8	none	none	100 mM Sodium acetate; pH 4.5	1 M di-Ammonium hydrogen phosphate
E9	none	none	100 mM MES; pH 6.5	1.6 M Magnesium sulfate
E10	10 % w/v Polyethylene glycol 6,000	none	100 mM BICINE; pH 9.0	none
E11	14.4 % w/v Polyethylene glycol 8,000	20 % v/v Glycerol	80 mM MES; pH 6.5	160 mM Calcium acetate
E12	10 % w/v Polyethylene glycol 8,000	none	100 mM Imidazole; pH 8.0	none
F1	30 % v/v Jeffamine® M-600	none	100 mM MES; pH 6.5	50 mM Cesium chloride
F2	none	none	100 mM tri-Sodium citrate; pH 5.0	3.15 M Ammonium sulfate
F3	20 % v/v 2-Methyl-2,4-pentanediol	none	100 mM TRIS; pH 8.0	none
F4	20 % v/v Jeffamine® M-600	none	100 mM HEPES; pH 7.5	none
F5	50 % v/v Ethylene glycol	none	100 mM TRIS; pH 8.5	200 mM Magnesium chloride
F6	10 % v/v 2-Methyl-2,4-pentanediol	none	100 mM BICINE; pH 9.0	none
F7	none	none	none	800 mM di-Sodium succinate; pH 7.0
F8	none	none	none	2.1 M di-Sodium DL-malate; pH 7.0
F9	none	none	none	2.4 M di-Sodium malonate; pH 7.0
F10	0.5 % w/v Jeffamine® ED-2001; pH 7.0	none	100 mM HEPES; pH 7.0	1.1 M di-Sodium malonate; pH 7.0
F11	1 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM HEPES; pH 7.0	1 M di-Sodium succinate; pH 7.0
F12	30 % v/v Jeffamine® M-600; pH 7.0	none	100 mM HEPES; pH 7.0	none

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

