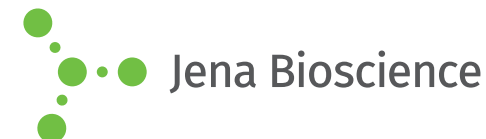




JBScreen JCSG++ HTS

Cat.-No.: CS-206L

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
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

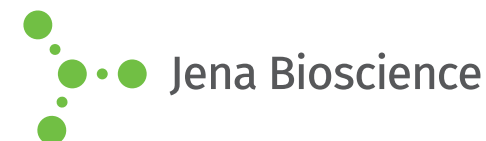




JBScreen JCSG++ HTS

Cat.-No.: CS-206L

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
C1	20 % w/v Polyethylene glycol 8,000	none	100 mM Potassium phosphate citrate; pH 4.2	200 mM Sodium chloride
C2	20 % w/v Polyethylene glycol 6,000	none	100 mM tri-Sodium citrate; pH 4.0	1 M Lithium chloride
C3	20 % w/v Polyethylene glycol 3,350	none	none	200 mM Ammonium nitrate; pH 6.3
C4	10 % w/v Polyethylene glycol 6,000	none	100 mM HEPES; pH 7.0	none
C5	800 mM Sodium di-hydrogen phosphate	800 mM Potassium di-hydrogen phosphate	100 mM HEPES; pH 7.5	none
C6	40 % v/v Polyethylene glycol 300	none	100 mM Potassium phosphate citrate; pH 4.2	none
C7	10 % w/v Polyethylene glycol 3,000	none	100 mM Sodium acetate; pH 4.5	200 mM Zinc acetate
C8	20 % v/v Ethanol	none	100 mM TRIS; pH 8.5	none
C9	25 % v/v 1,2-Propanediol	10 % v/v Glycerol	100 mM Sodium Potassium phosphate; pH 6.2	none
C10	10 % w/v Polyethylene glycol 20,000	2 % v/v 1,4-Dioxane	100 mM BICINE; pH 9.0	none
C11	none	none	100 mM Sodium acetate; pH 4.6	2 M Ammonium sulfate
C12	10 % w/v Polyethylene glycol 1,000	10 % w/v Polyethylene glycol 8,000	none	none
D1	24 % w/v Polyethylene glycol 1,500	20 % v/v Glycerol	none	none
D2	30 % v/v Polyethylene glycol 400	none	100 mM HEPES; pH 7.5	200 mM Magnesium chloride
D3	50 % v/v Polyethylene glycol 200	none	100 mM Sodium Potassium phosphate; pH 6.2	200 mM Sodium chloride
D4	30 % w/v Polyethylene glycol 8,000	none	100 mM Sodium acetate; pH 4.5	200 mM Lithium sulfate
D5	70 % v/v 2-Methyl-2,4-pentanediol	none	100 mM HEPES; pH 7.5	none
D6	20 % w/v Polyethylene glycol 8,000	none	100 mM TRIS; pH 8.5	200 mM Magnesium chloride
D7	40 % v/v Polyethylene glycol 400	none	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
D8	40 % v/v 2-Methyl-2,4-pentanediol	none	100 mM TRIS; pH 8.0	none
D9	25.5 % w/v Polyethylene glycol 4,000	15 % v/v Glycerol	none	170 mM Ammonium sulfate
D10	40 % v/v Polyethylene glycol 300	none	100 mM MES; pH 6.5	200 mM Calcium acetate
D11	14 % v/v 2-Propanol	30 % v/v Glycerol	70 mM Sodium acetate; pH 4.6	140 mM Calcium chloride
D12	16 % w/v Polyethylene glycol 8,000	20 % v/v Glycerol	none	40 mM Potassium di-hydrogen phosphate

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





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



No.	Precipitant	Precipitant 2	Buffer	Additive
G1	30 % w/v Jeffamine® ED-2001; pH 7.0	none	100 mM HEPES; pH 7.0	none
G2	22 % w/v Poly(acrylic acid sodium salt) 5,100	none	100 mM HEPES; pH 7.5	20 mM Magnesium chloride
G3	20 % w/v Polyvinylpyrrolidone K15	none	100 mM TRIS; pH 8.5	100 mM Cobalt (II) chloride
G4	20 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM TRIS; pH 8.5	200 mM Trimethylamine N-oxide
G5	12 % w/v Polyethylene glycol 3,350	none	100 mM HEPES; pH 7.5	5 mM Cobalt (II) chloride, 5 mM Cadmium chloride, 5 mM Nickel (II) chloride, 5 mM Magnesium chloride
G6	20 % w/v Polyethylene glycol 3,350	none	none	200 mM di-Sodium malonate; pH 7.0
G7	15 % w/v Polyethylene glycol 3,350	none	none	100 mM di-Sodium succinate; pH 7.0
G8	20 % w/v Polyethylene glycol 3,350	none	none	150 mM di-Sodium DL-malate; pH 7.0
G9	30 % w/v Polyethylene glycol monomethyl ether 2,000	none	none	100 mM Potassium thiocyanate
G10	30 % w/v Polyethylene glycol monomethyl ether 2,000	none	none	150 mM Potassium bromide
G11	none	none	100 mM BIS-TRIS; pH 5.5	2 M Ammonium sulfate
G12	none	none	100 mM BIS-TRIS; pH 5.5	3 M Sodium chloride
H1	none	none	100 mM BIS-TRIS; pH 5.5	300 mM Magnesium formate
H2	1 % w/v Polyethylene glycol 3,350	none	100 mM BIS-TRIS; pH 5.5	1 M Ammonium sulfate
H3	25 % w/v Polyethylene glycol 3,350	none	100 mM BIS-TRIS; pH 5.5	none
H4	45 % v/v 2-Methyl-2,4-pentanediol	none	100 mM BIS-TRIS; pH 5.5	200 mM Calcium chloride
H5	45 % v/v 2-Methyl-2,4-pentanediol	none	100 mM BIS-TRIS; pH 5.5	200 mM Ammonium acetate
H6	17 % w/v Polyethylene glycol 10,000	none	100 mM BIS-TRIS; pH 5.5	100 mM Ammonium acetate
H7	25 % w/v Polyethylene glycol 3,350	none	100 mM BIS-TRIS; pH 5.5	200 mM Ammonium sulfate
H8	25 % w/v Polyethylene glycol 3,350	none	100 mM BIS-TRIS; pH 5.5	200 mM Sodium chloride
H9	25 % w/v Polyethylene glycol 3,350	none	100 mM BIS-TRIS; pH 5.5	200 mM Lithium sulfate
H10	25 % w/v Polyethylene glycol 3,350	none	100 mM BIS-TRIS; pH 5.5	200 mM Ammonium acetate
H11	25 % w/v Polyethylene glycol 3,350	none	100 mM BIS-TRIS; pH 5.5	200 mM Magnesium chloride
H12	45 % v/v 2-Methyl-2,4-pentanediol	none	100 mM HEPES; pH 7.5	200 mM Ammonium acetate

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