



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



## JBScreen Basic HTS

Cat.-No.: CS-203L

# SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
C1	8 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	none
C2	20 % w/v Polyethylene glycol 4,000	20 % v/v 2-Propanol	100 mM tri-Sodium citrate; pH 5.6	none
C3	20 % w/v Polyethylene glycol 4,000	10 % v/v 2-Propanol	100 mM HEPES; pH 7.5	none
C4	25 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate
C5	30 % w/v Polyethylene glycol 4,000	none	none	200 mM Ammonium sulfate
C6	30 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	200 mM Ammonium acetate
C7	30 % w/v Polyethylene glycol 4,000	none	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium acetate
C8	30 % w/v Polyethylene glycol 4,000	none	100 mM TRIS; pH 8.5	200 mM Sodium acetate
C9	30 % w/v Polyethylene glycol 4,000	none	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
C10	30 % w/v Polyethylene glycol 4,000	none	100 mM TRIS; pH 8.5	200 mM Magnesium chloride
C11	30 % w/v Polyethylene glycol monomethyl ether 5,000	none	100 mM MES; pH 6.5	200 mM Ammonium sulfate
C12	10 % w/v Polyethylene glycol 6,000	2 M Sodium chloride	none	none
D1	10 % w/v Polyethylene glycol 6,000	5 % v/v 2-Methyl-2,4-pentanediol	100 mM HEPES; pH 7.5	none
D2	2 % w/v Polyethylene glycol 8,000	1 M Lithium sulfate	none	none
D3	8 % w/v Polyethylene glycol 8,000	none	100 mM TRIS; pH 8.5	none
D4	10 % w/v Polyethylene glycol 8,000	8 % v/v Ethylene glycol	100 mM HEPES; pH 7.5	none
D5	15 % w/v Polyethylene glycol 8,000	500 mM Lithium sulfate	none	none
D6	18 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Calcium acetate
D7	18 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Zinc acetate
D8	20 % w/v Polyethylene glycol 8,000	none	none	50 mM Potassium di-hydrogen phosphate
D9	20 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Magnesium acetate
D10	30 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Sodium acetate
D11	30 % w/v Polyethylene glycol 8,000	none	none	200 mM Ammonium sulfate
D12	30 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Ammonium sulfate

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

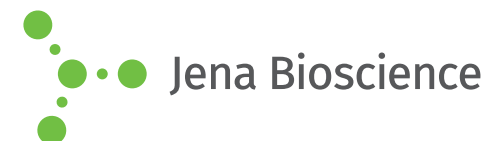




## JBScreen Basic HTS

Cat.-No.: CS-203L

# SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
E1	10 % w/v Polyethylene glycol 10,000	2 % v/v 1,4-Dioxane	100 mM BICINE; pH 9.5	none
E2	20 % w/v Polyethylene glycol 10,000	none	100 mM HEPES; pH 7.5	none
E3	12 % w/v Polyethylene glycol 20,000	none	100 mM MES; pH 6.5	none
E4	5 % v/v 2-Propanol	2 M Ammonium sulfate	none	none
E5	20 % v/v 2-Propanol	none	100 mM HEPES; pH 7.5	200 mM tri-Sodium citrate
E6	20 % v/v 2-Propanol	none	100 mM Sodium acetate; pH 4.6	200 mM Calcium chloride
E7	30 % v/v 2-Propanol	none	100 mM HEPES; pH 7.5	200 mM Magnesium chloride
E8	30 % v/v 2-Propanol	none	100 mM TRIS; pH 8.5	200 mM Ammonium acetate
E9	10 % v/v 1,4-Dioxane	1.6 M Ammonium sulfate	100 mM MES; pH 6.5	none
E10	35 % v/v 1,4-Dioxane	none	none	none
E11	10 % v/v Ethanol	1.5 M Sodium chloride	none	none
E12	20 % v/v Ethanol	none	100 mM TRIS; pH 8.5	none
F1	25 % v/v 2-Methyl-2-propanol	none	100 mM TRIS; pH 8.5	none
F2	35 % v/v 2-Methyl-2-propanol	none	100 mM tri-Sodium citrate; pH 5.6	none
F3	1 M Imidazole; pH 7.0	none	none	none
F4	1 M Lithium sulfate	none	100 mM TRIS; pH 8.5	10 mM Nickel (II) chloride
F5	1.5 M Lithium sulfate	none	100 mM HEPES; pH 7.5	none
F6	400 mM Potassium Sodium tartrate	none	none	none
F7	800 mM Potassium Sodium tartrate	none	100 mM HEPES; pH 7.5	none
F8	1.4 M tri-Sodium citrate	none	100 mM HEPES; pH 7.5	none
F9	1.6 M tri-Sodium citrate; pH 6.5	none	none	none
F10	10 % v/v Jeffamine® M-600	none	100 mM tri-Sodium citrate; pH 5.6	10 mM Iron (III) chloride
F11	20 % v/v Jeffamine® M-600	none	100 mM HEPES; pH 7.5	none
F12	30 % v/v Jeffamine® M-600	none	100 mM MES; pH 6.5	50 mM Cesium chloride

\*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	800 mM Potassium di-hydrogen phosphate	800 mM Sodium di-hydrogen phosphate	100 mM HEPES; pH 7.5	none
G2	400 mM Ammonium di-hydrogen phosphate	none	none	none
G3	1 M Ammonium di-hydrogen phosphate	none	100 mM tri-Sodium citrate; pH 5.6	none
G4	2 M Ammonium di-hydrogen phosphate	none	100 mM TRIS; pH 8.5	none
G5	2 M Ammonium formate	none	100 mM Sodium acetate; pH 4.6	none
G6	4 M Ammonium formate	none	100 mM HEPES; pH 7.5	none
G7	2 M Ammonium formate	none	none	none
G8	500 mM Ammonium sulfate	1 M Lithium sulfate	100 mM tri-Sodium citrate; pH 5.6	none
G9	1.6 M Ammonium sulfate	none	100 mM HEPES; pH 7.5	100 mM Sodium chloride
G10	1.8 M Ammonium sulfate	none	100 mM MES; pH 6.5	10 mM Cobalt (II) chloride
G11	2 M Ammonium sulfate	none	100 mM TRIS; pH 8.5	none
G12	2 M Ammonium sulfate	none	none	none
H1	2 M Ammonium sulfate	none	100 mM Sodium acetate; pH 4.6	none
H2	2 M Ammonium sulfate	none	100 mM tri-Sodium citrate; pH 5.6	200 mM Potassium Sodium tartrate
H3	200 mM Magnesium formate	none	none	none
H4	1.6 M Magnesium sulfate	none	100 mM MES; pH 6.5	none
H5	2 M Magnesium chloride	none	100 mM BICINE; pH 9.5	none
H6	1 M Sodium acetate	none	100 mM Imidazole; pH 6.5	none
H7	1 M Sodium acetate	none	100 mM HEPES; pH 7.5	50 mM Cadmium sulfate
H8	1.4 M Sodium acetate	none	100 mM MES; pH 6.5	none
H9	500 mM Sodium chloride	10 mM Magnesium chloride	none	10 mM Cetyltrimethylammonium bromide
H10	2 M Sodium chloride	none	100 mM Sodium acetate; pH 4.6	none
H11	2 M Sodium chloride	none	100 mM MES; pH 6.5	100 mM Sodium di-hydrogen phosphate, 100 mM Potassium di-hydrogen phosphate
H12	4.3 M Sodium chloride	none	100 mM HEPES; pH 7.5	none

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