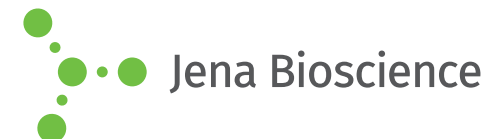




JBScreen XP Up

Kit including JBScreen XP and TEW
Cat.-No.: CS-351

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	50 mM 1,6-Hexanediol	none	100 mM Sodium acetate; pH 4.6	5 mM Cobalt (II) chloride
A4	1 M 1,6-Hexanediol	none	100 mM tri-Sodium citrate; pH 5.6	none
A5	1 M 1,6-Hexanediol	none	100 mM HEPES; pH 7.5	10 mM Magnesium chloride
A6	5 % v/v 2-Methyl-2,4-pentanediol	none	100 mM Sodium acetate; pH 4.6	2 mM Calcium chloride
A7	5 % v/v 2-Methyl-2,4-pentanediol	none	100 mM Sodium acetate; pH 4.6	10 mM Sodium chloride
A8	10 % v/v 2-Methyl-2,4-pentanediol	none	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium acetate
A9	10 % v/v 2-Methyl-2,4-pentanediol	none	100 mM HEPES; pH 7.5	10 mM Ammonium sulfate
A10	10 % v/v 2-Methyl-2,4-pentanediol	none	100 mM TRIS; pH 8.5	200 mM Ammonium di-hydrogen phosphate
A11	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM HEPES; pH 7.5	none
A12	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM HEPES; pH 7.5	100 mM tri-Sodium citrate
B1	2 % v/v Polyethylene glycol 400	none	100 mM tri-Sodium citrate; pH 5.6	none
B2	2 % v/v Polyethylene glycol 400	none	100 mM tri-Sodium citrate; pH 5.6	500 mM Sodium chloride
B3	2 % v/v Polyethylene glycol 400	none	100 mM HEPES; pH 7.5	10 mM Calcium chloride
B4	2 % v/v Polyethylene glycol 400	1 M Ammonium sulfate	100 mM HEPES; pH 7.5	none
B5	5 % v/v Polyethylene glycol 400	none	100 mM Sodium acetate; pH 4.6	10 mM Calcium chloride
B6	30 % v/v Polyethylene glycol 400	none	100 mM HEPES; pH 7.5	200 mM Magnesium chloride
B7	10 % v/v Polyethylene glycol monomethyl ether 550	none	100 mM MES; pH 6.5	none
B8	20 % v/v Polyethylene glycol monomethyl ether 550	none	100 mM HEPES; pH 7.5	100 mM Sodium chloride
B9	10 % w/v Polyethylene glycol 1,000	5 % w/v Polyethylene glycol 8,000	100 mM tri-Sodium citrate; pH 5.6	none
B10	30 % w/v Polyethylene glycol 1,500	none	100 mM HEPES; pH 7.5	none
B11	5 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM TRIS; pH 8.5	none
B12	15 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM Sodium acetate; pH 4.6	10 mM Ammonium sulfate
C1	20 % w/v Polyethylene glycol monomethyl ether 2,000	none	none	none

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

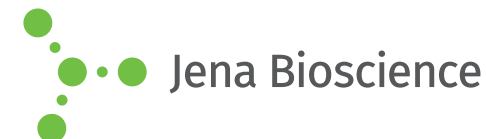




JBScreen XP Up

Kit including JBScreen XP and TEW
Cat.-No.: CS-351

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
C2	8 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	none
C3	10 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	10 mM Ammonium sulfate
C4	10 % w/v Polyethylene glycol 4,000	5 % v/v 2-Propanol	100 mM tri-Sodium citrate; pH 5.6	none
C5	15 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	10 mM Ammonium acetate
C6	20 % w/v Polyethylene glycol 4,000	none	100 mM TRIS; pH 8.5	10 mM Magnesium chloride
C7	20 % w/v Polyethylene glycol 4,000	none	100 mM TRIS; pH 8.5	200 mM Sodium acetate
C8	20 % w/v Polyethylene glycol 4,000	5 % v/v 2-Propanol	100 mM HEPES; pH 7.5	none
C9	30 % w/v Polyethylene glycol 4,000	none	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium acetate
C10	30 % w/v Polyethylene glycol 4,000	none	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
C11	10 % w/v Polyethylene glycol monomethyl ether 5,000	none	100 mM MES; pH 6.5	10 mM Ammonium sulfate
C12	10 % w/v Polyethylene glycol 6,000	none	100 mM tri-Sodium citrate; pH 5.6	100 mM Sodium chloride
D1	10 % w/v Polyethylene glycol 6,000	5 % v/v 2-Methyl-2,4-pentanediol	100 mM HEPES; pH 7.5	none
D2	5 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	2 mM Calcium acetate
D3	8 % w/v Polyethylene glycol 8,000	none	100 mM TRIS; pH 8.5	none
D4	10 % w/v Polyethylene glycol 8,000	none	none	10 mM Ammonium sulfate
D5	10 % w/v Polyethylene glycol 8,000	none	none	10 mM Potassium di-hydrogen phosphate
D6	10 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	5 mM Ammonium sulfate
D7	10 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	5 mM Sodium acetate
D8	10 % w/v Polyethylene glycol 8,000	8 % v/v Ethylene glycol	100 mM HEPES; pH 7.5	none
D9	15 % w/v Polyethylene glycol 8,000	500 mM Lithium sulfate	100 mM Sodium acetate; pH 4.6	none
D10	18 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	none
D11	20 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Magnesium acetate
D12	10 % w/v Polyethylene glycol 10,000	2 % v/v 1,4-Dioxane	100 mM tri-Sodium citrate; pH 5.6	none
E1	20 % w/v Polyethylene glycol 10,000	none	100 mM HEPES; pH 7.5	none
E2	12 % w/v Polyethylene glycol 20,000	none	100 mM MES; pH 6.5	none

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

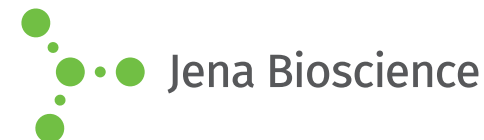




JBScreen XP Up

Kit including JBScreen XP and TEW
Cat.-No.: CS-351

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
E3	5 % v/v 2-Propanol	1 M Ammonium sulfate	none	none
E4	10 % v/v 2-Propanol	none	100 mM Sodium acetate; pH 4.6	none
E5	10 % v/v 2-Propanol	none	100 mM TRIS; pH 8.5	10 mM Ammonium acetate
E6	20 % v/v 2-Propanol	none	100 mM HEPES; pH 7.5	75 mM Magnesium chloride
E7	20 % v/v 2-Propanol	none	100 mM HEPES; pH 7.5	100 mM tri-Sodium citrate
E8	10 % v/v 1,4-Dioxane	none	100 mM MES; pH 6.5	20 mM Ammonium sulfate
E9	25 % v/v 1,4-Dioxane	none	none	none
E10	10 % v/v Ethanol	none	none	100 mM Sodium chloride
E11	20 % v/v Ethanol	none	100 mM TRIS; pH 8.5	none
E12	10 % v/v 2-Methyl-2-propanol	none	100 mM HEPES; pH 7.5	none
F1	25 % v/v 2-Methyl-2-propanol	none	100 mM TRIS; pH 8.5	none
F2	500 mM Imidazole; pH 7.0	none	none	none
F3	100 mM Lithium sulfate	none	100 mM TRIS; pH 8.5	none
F4	1.5 M Lithium sulfate	none	100 mM HEPES; pH 7.5	none
F5	400 mM Potassium Sodium tartrate	none	none	none
F6	400 mM Potassium Sodium tartrate	none	100 mM Imidazole; pH 6.5	none
F7	400 mM Potassium Sodium tartrate	none	100 mM BICINE; pH 9.5	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	5 % v/v Jeffamine® M-600; pH 7.0	none	100 mM tri-Sodium citrate; pH 5.6	10 mM Iron (III) chloride
F11	20 % v/v Jeffamine® M-600; pH 7.0	none	100 mM MES; pH 6.5	10 mM Cesium chloride
F12	20 % v/v Jeffamine® M-600; pH 7.0	none	100 mM HEPES; pH 7.5	none
G1	100 mM Potassium di-hydrogen phosphate	none	100 mM tri-Sodium citrate; pH 5.6	100 mM Sodium di-hydrogen phosphate
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

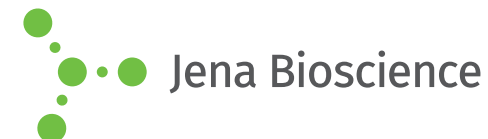
*pH values indicated are those of the 1.0 M buffer stock solution prior to dilution with other components



JBScreen XP Up

Kit including JBScreen XP and TEW
Cat.-No.: CS-351

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
G4	2 M Ammonium di-hydrogen phosphate	none	100 mM TRIS; pH 8.5	none
G5	1.5 M Ammonium formate	none	100 mM Sodium acetate; pH 4.6	none
G6	2 M Ammonium formate	none	none	none
G7	4 M Ammonium formate	none	100 mM HEPES; pH 7.5	none
G8	500 mM Ammonium sulfate	1 M Lithium sulfate	100 mM tri-Sodium citrate; pH 5.6	none
G9	1 M Ammonium sulfate	none	100 mM MES; pH 6.5	10 mM Cobalt (II) chloride
G10	1 M Ammonium sulfate	none	100 mM HEPES; pH 7.5	100 mM Sodium chloride
G11	2 M Ammonium sulfate	none	none	none
G12	2 M Ammonium sulfate	none	100 mM Sodium acetate; pH 4.6	none
H1	2 M Ammonium sulfate	none	100 mM tri-Sodium citrate; pH 5.6	200 mM Potassium Sodium tartrate
H2	2 M Ammonium sulfate	none	100 mM TRIS; pH 8.5	none
H3	200 mM Magnesium formate	none	none	none
H4	1.6 M Magnesium sulfate	none	100 mM MES; pH 6.5	none
H5	1.2 M Magnesium chloride	none	100 mM BICINE; pH 9.5	none
H6	100 mM Sodium acetate	none	100 mM tri-Sodium citrate; pH 5.6	2 mM Cadmium sulfate
H7	500 mM Sodium acetate	none	100 mM MES; pH 6.5	none
H8	1 M Sodium acetate	none	100 mM Imidazole; pH 6.5	none
H9	100 mM Sodium chloride	none	none	10 mM Magnesium chloride
H10	1 M Sodium chloride	none	100 mM Sodium acetate; pH 4.6	none
H11	1 M Sodium chloride	none	100 mM HEPES; pH 7.5	none
H12	2 M Sodium chloride	none	100 mM MES; pH 6.5	100 mM Sodium di-hydrogen phosphate, 100 mM Potassium di-hydrogen phosphate

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

