

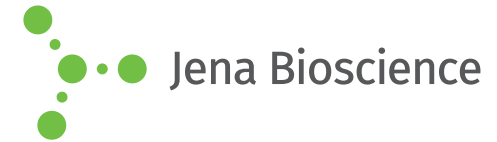


JBScreen Classic HTS I

(PEG based)

Cat.-No.: CS-201L

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive	Classic Bulk No.
A1	15.00 % w/v Polyethylene glycol 400	None	100 mM Sodium acetate; pH 4.6	100 mM Calcium chloride	1/A1
A2	15.00 % w/v Polyethylene glycol 400	None	100 mM HEPES; pH 7.5	200 mM Magnesium chloride	1/A3
A3	25.00 % w/v Polyethylene glycol 400	None	100 mM Sodium acetate; pH 4.6	100 mM Magnesium chloride	1/A5
A4	25.00 % w/v Polyethylene glycol 400	None	100 mM TRIS; pH 8.5	200 mM Lithium sulfate	1/A6
A5	30.00 % w/v Polyethylene glycol 400	None	100 mM Sodium acetate; pH 4.6	100 mM Calcium chloride	1/B2
A6	30.00 % w/v Polyethylene glycol 400	None	100 mM MES; pH 6.5	100 mM Sodium acetate	1/B3
A7	30.00 % w/v Polyethylene glycol 400	None	100 mM HEPES; pH 7.5	200 mM Magnesium chloride	1/B5
A8	30.00 % w/v Polyethylene glycol 400	None	100 mM TRIS; pH 8.5	200 mM tri-Sodium citrate	1/B6
A9	30.00 % w/v Polyethylene glycol monomethyl ether 550	None	100 mM BICINE; pH 9.0	100 mM Sodium chloride	1/C1
A10	25.00 % w/v Polyethylene glycol monomethyl ether 550	None	100 mM MES; pH 6.5	10.00 mM Zinc sulfate	1/C2
A11	25.00 % w/v Polyethylene glycol 1,000	None	100 mM HEPES; pH 7.5	None	1/C3
A12	30.00 % w/v Polyethylene glycol 1,000	None	100 mM TRIS; pH 8.5	None	1/C4
B1	15.00 % w/v Polyethylene glycol 1,500	None	None	None	1/C5
B2	20.00 % w/v Polyethylene glycol 1,500	None	100 mM HEPES; pH 7.5	None	1/C6
B3	30.00 % w/v Polyethylene glycol 1,500	None	None	None	1/D1
B4	20.00 % w/v Polyethylene glycol monomethyl ether 2,000	None	100 mM TRIS; pH 8.5	10.00 mM Nickel (II) chloride	1/D2
B5	25.00 % w/v Polyethylene glycol monomethyl ether 2,000	None	None	None	1/D3
B6	20.00 % w/v Polyethylene glycol 3,000	None	100 mM HEPES; pH 7.5	200 mM Sodium acetate	1/D5
B7	30.00 % w/v Polyethylene glycol 3,000	None	100 mM TRIS; pH 8.5	200 mM Lithium sulfate	1/D6
B8	4.00 % w/v Polyethylene glycol 4,000	None	100 mM Sodium acetate; pH 4.6	None	2/A1
B9	8.00 % w/v Polyethylene glycol 4,000	None	None	None	2/A2
B10	8.00 % w/v Polyethylene glycol 4,000	None	100 mM Sodium acetate; pH 4.6	None	2/A3
B11	10.00 % w/v Polyethylene glycol 4,000	None	100 mM MES; pH 6.5	200 mM Magnesium chloride	2/A4
B12	12.00 % w/v Polyethylene glycol 4,000	None	100 mM HEPES; pH 7.5	100 mM Sodium acetate	2/A5

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



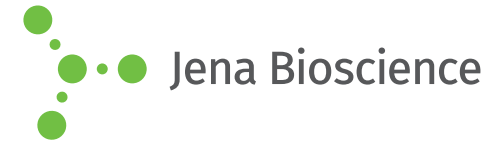


JBScreen Classic HTS I

(PEG based)

Cat.-No.: CS-201L

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive	Classic Bulk No.
C1	16.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	200 mM Lithium sulfate	2/B1
C2	16.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	200 mM Sodium acetate	2/B2
C3	18.00 % w/v Polyethylene glycol 4,000	None	100 mM Sodium acetate; pH 4.6	None	2/B4
C4	20.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	200 mM Lithium sulfate	2/B5
C5	20.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	200 mM Calcium chloride	2/B6
C6	25.00 % w/v Polyethylene glycol 4,000	None	100 mM Sodium acetate; pH 4.6	None	2/C2
C7	25.00 % w/v Polyethylene glycol 4,000	None	100 mM MES; pH 6.5	200 mM Magnesium chloride	2/C3
C8	25.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	200 mM Calcium chloride	2/C4
C9	30.00 % w/v Polyethylene glycol 4,000	None	None	None	2/C5
C10	30.00 % w/v Polyethylene glycol 4,000	None	100 mM Sodium acetate; pH 4.6	100 mM Magnesium chloride	2/C6
C11	30.00 % w/v Polyethylene glycol 4,000	None	100 mM HEPES; pH 7.5	200 mM Calcium chloride	2/D2
C12	30.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	200 mM Sodium acetate	2/D4
D1	30.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	200 mM Magnesium chloride	2/D5
D2	35.00 % w/v Polyethylene glycol 4,000	None	None	None	2/D6
D3	8.00 % w/v Polyethylene glycol 4,000	800 mM Lithium chloride	100 mM TRIS; pH 8.5	None	3/A1
D4	10.00 % w/v Polyethylene glycol 4,000	20.00 % w/v 2-Propanol	None	None	3/A2
D5	10.00 % w/v Polyethylene glycol 4,000	10.00 % w/v 2-Propanol	100 mM tri-Sodium citrate; pH 5.6	None	3/A3
D6	10.00 % w/v Polyethylene glycol 4,000	20.00 % w/v 2-Propanol	100 mM HEPES; pH 7.5	None	3/A5
D7	12.00 % w/v Polyethylene glycol 4,000	None	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate	3/A6
D8	15.00 % w/v Polyethylene glycol 4,000	None	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium sulfate	3/B2
D9	16.00 % w/v Polyethylene glycol 4,000	10.00 % w/v 2-Propanol	100 mM HEPES; pH 7.5	200 mM Ammonium sulfate	3/B3
D10	20.00 % w/v Polyethylene glycol 4,000	None	None	200 mM Ammonium sulfate	3/B4
D11	20.00 % w/v Polyethylene glycol 4,000	10.00 % w/v Glycerol	None	200 mM Magnesium sulfate	3/B5
D12	20.00 % w/v Polyethylene glycol 4,000	20.00 % w/v 2-Propanol	None	100 mM tri-Sodium citrate	3/C1

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

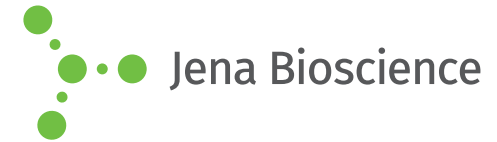


JBScreen Classic HTS I

(PEG based)

Cat.-No.: CS-201L

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive	Classic Bulk No.
E1	20.00 % w/v Polyethylene glycol 4,000	None	100 mM MES; pH 6.5	600 mM Sodium chloride	3/C2
E2	20.00 % w/v Polyethylene glycol 4,000	10.00 % w/v 2-Propanol	100 mM HEPES; pH 7.5	None	3/C3
E3	22.00 % w/v Polyethylene glycol 4,000	None	None	100 mM Sodium acetate, 200 mM Ammonium sulfate	3/C4
E4	25.00 % w/v Polyethylene glycol 4,000	None	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium sulfate	3/C6
E5	25.00 % w/v Polyethylene glycol 4,000	200 mM Lithium sulfate	100 mM HEPES; pH 7.5	100 mM Sodium acetate	3/D1
E6	25.00 % w/v Polyethylene glycol 4,000	8.00 % w/v 2-Propanol	None	100 mM Sodium acetate	3/D2
E7	30.00 % w/v Polyethylene glycol 4,000	None	None	200 mM Ammonium sulfate	3/D3
E8	30.00 % w/v Polyethylene glycol 4,000	None	100 mM tri-Sodium citrate; pH 5.6	100 mM Ammonium sulfate	3/D5
E9	32.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	800 mM Lithium chloride	3/D6
E10	25.00 % w/v Polyethylene glycol monomethyl ether 5,000	None	100 mM TRIS; pH 8.5	200 mM Lithium sulfate	4/A1
E11	30.00 % w/v Polyethylene glycol monomethyl ether 5,000	None	100 mM MES; pH 6.5	200 mM Ammonium sulfate	4/A2
E12	3.00 % w/v Polyethylene glycol 6,000	None	100 mM TRIS; pH 8.5	100 mM Potassium chloride	4/A3
F1	10.00 % w/v Polyethylene glycol 6,000	None	None	10.00 mM Magnesium chloride	4/A4
F2	12.00 % w/v Polyethylene glycol 6,000	2.00 M Sodium chloride	None	None	4/A5
F3	15.00 % w/v Polyethylene glycol 6,000	5.00 % w/v Glycerol	None	None	4/A6
F4	15.00 % w/v Polyethylene glycol 6,000	50.00 mM Potassium chloride	None	10.00 mM Magnesium chloride	4/B1
F5	20.00 % w/v Polyethylene glycol 6,000	None	50.00 mM Imidazole; pH 8.0	None	4/B3
F6	25.00 % w/v Polyethylene glycol 6,000	None	100 mM HEPES; pH 7.5	100 mM Lithium chloride	4/B4
F7	28.00 % w/v Polyethylene glycol 6,000	500 mM Lithium chloride	100 mM TRIS; pH 8.5	None	4/B5
F8	30.00 % w/v Polyethylene glycol 6,000	1.00 M Lithium chloride	None	100 mM Sodium acetate	4/B6
F9	2.00 % w/v Polyethylene glycol 8,000	500 mM Lithium sulfate	None	None	4/C2
F10	2.00 % w/v Polyethylene glycol 8,000	1.00 M Lithium sulfate	None	None	4/C3
F11	4.00 % w/v Polyethylene glycol 8,000	None	None	None	4/C4
F12	8.00 % w/v Polyethylene glycol 8,000	200 mM Lithium chloride	None	50.00 mM Magnesium sulfate	4/C5

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



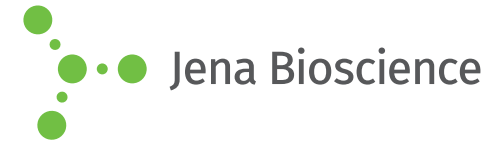


JBScreen Classic HTS I

(PEG based)

Cat.-No.: CS-201L

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive	Classic Bulk No.
G1	10.00 % w/v Polyethylene glycol 8,000	None	100 mM MES; pH 6.5	200 mM Zinc acetate	4/D1
G2	10.00 % w/v Polyethylene glycol 8,000	None	100 mM HEPES; pH 7.5	200 mM Calcium acetate	4/D2
G3	10.00 % w/v Polyethylene glycol 8,000	None	None	50.00 mM Magnesium acetate, 100 mM Sodium acetate	4/D3
G4	10.00 % w/v Polyethylene glycol 8,000	None	None	200 mM Magnesium acetate	4/D4
G5	10.00 % w/v Polyethylene glycol 8,000	10.00 % w/v Ethylene glycol	100 mM HEPES; pH 7.5	None	4/D5
G6	12.00 % w/v Polyethylene glycol 8,000	10.00 % w/v Glycerol	None	500 mM Potassium chloride	5/A2
G7	15.00 % w/v Polyethylene glycol 8,000	None	None	200 mM Ammonium sulfate	5/A3
G8	15.00 % w/v Polyethylene glycol 8,000	500 mM Lithium sulfate	None	None	5/A4
G9	15.00 % w/v Polyethylene glycol 8,000	None	100 mM MES; pH 6.5	200 mM Sodium acetate	5/A5
G10	18.00 % w/v Polyethylene glycol 8,000	None	100 mM HEPES; pH 7.5	200 mM Calcium acetate	5/B1
G11	18.00 % w/v Polyethylene glycol 8,000	2.00 % w/v 2-Propanol	100 mM HEPES; pH 7.5	100 mM Sodium acetate	5/B2
G12	18.00 % w/v Polyethylene glycol 8,000	None	100 mM TRIS; pH 8.5	200 mM Lithium sulfate	5/B3
H1	20.00 % w/v Polyethylene glycol 8,000	None	100 mM MES; pH 6.5	200 mM Magnesium acetate	5/B5
H2	20.00 % w/v Polyethylene glycol 8,000	None	100 mM CHES; pH 9.5	None	5/B6
H3	25.00 % w/v Polyethylene glycol 8,000	None	None	200 mM Lithium chloride	5/C2
H4	30.00 % w/v Polyethylene glycol 8,000	None	None	200 mM Ammonium sulfate	5/C3
H5	8.00 % w/v Polyethylene glycol 10,000	None	100 mM Sodium acetate; pH 4.6	None	5/C4
H6	14.00 % w/v Polyethylene glycol 10,000	None	100 mM Imidazole; pH 8.0	None	5/C5
H7	18.00 % w/v Polyethylene glycol 10,000	20.00 % w/v Glycerol	100 mM TRIS; pH 8.5	100 mM Sodium chloride	5/D1
H8	20.00 % w/v Polyethylene glycol 10,000	None	100 mM HEPES; pH 7.5	None	5/D2
H9	30.00 % w/v Polyethylene glycol 10,000	None	100 mM TRIS; pH 8.5	None	5/D3
H10	10.00 % w/v Polyethylene glycol 20,000	None	100 mM MES; pH 6.5	None	5/D4
H11	17.00 % w/v Polyethylene glycol 20,000	None	100 mM TRIS; pH 8.5	100 mM Magnesium chloride	5/D5
H12	20.00 % w/v Polyethylene glycol 20,000	None	None	None	5/D6

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

