

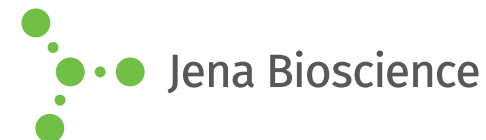


JBScreen Classic HTS II

(Ammonium Sulfate, MPD, Alcohol and Salt based)

Cat.-No.: CS-202L

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive	Classic Bulk No.
A1	500 mM Ammonium sulfate	1.00 M Lithium sulfate	None	100 mM tri-Sodium citrate	6/A1
A2	1.00 M Ammonium sulfate	None	100 mM Sodium acetate; pH 4.6	None	6/A3
A3	1.00 M Ammonium sulfate	2.00 % w/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	None	6/A4
A4	1.00 M Ammonium sulfate	None	100 mM TRIS; pH 8.5	None	6/A5
A5	1.2 M Ammonium sulfate	3.00 % w/v 2-Propanol	None	50.00 mM tri-Sodium citrate	6/A6
A6	1.5 M Ammonium sulfate	15.00 % w/v Glycerol	100 mM TRIS; pH 8.5	None	6/B1
A7	1.6 M Ammonium sulfate	1.00 M Lithium sulfate	None	None	6/B3
A8	1.6 M Ammonium sulfate	None	100 mM HEPES; pH 7.5	200 mM Sodium chloride	6/B4
A9	1.6 M Ammonium sulfate	2.00 % w/v Polyethylene glycol 1,000	100 mM HEPES; pH 7.5	None	6/B5
A10	1.8 M Ammonium sulfate	None	100 mM MES; pH 6.5	None	6/B6
A11	2.00 M Ammonium sulfate	2.00 M Sodium chloride	None	None	6/C1
A12	2.00 M Ammonium sulfate	None	100 mM Sodium acetate; pH 4.6	None	6/C2
B1	2.00 M Ammonium sulfate	5.00 % w/v Polyethylene glycol 400	100 mM MES; pH 6.5	None	6/C3
B2	2.00 M Ammonium sulfate	None	100 mM TRIS; pH 8.5	None	6/C4
B3	2.2 M Ammonium sulfate	20.00 % w/v Glycerol	None	None	6/C6
B4	2.4 M Ammonium sulfate	None	None	100 mM tri-Sodium citrate	6/D1
B5	3.00 M Ammonium sulfate	1.00 % w/v 2-Methyl-2,4-pentanediol	None	None	6/D2
B6	3.00 M Ammonium sulfate	10.00 % w/v Glycerol	None	None	6/D3
B7	3.5 M Ammonium sulfate	None	100 mM HEPES; pH 7.5	None	6/D4
B8	3.5 M Ammonium sulfate	1.00 % w/v 2-Methyl-2,4-pentanediol	100 mM MES; pH 6.5	None	6/D5
B9	10.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM HEPES; pH 7.5	100 mM tri-Sodium citrate	7/A1
B10	12.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM TRIS; pH 8.5	50.00 mM Magnesium chloride	7/A2
B11	15.00 % w/v 2-Methyl-2,4-pentanediol	5.00 % w/v Polyethylene glycol 4,000	100 mM Imidazole; pH 8.0	None	7/A4
B12	15.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium acetate	7/A5

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



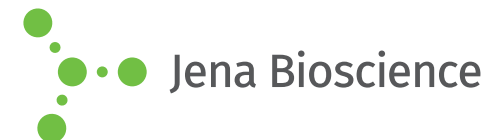


JBScreen Classic HTS II

(Ammonium Sulfate, MPD, Alcohol and Salt based)

Cat.-No.: CS-202L

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive	Classic Bulk No.
C1	15.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM HEPES; pH 7.5	200 mM tri-Sodium citrate	7/B1
C2	20.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM HEPES; pH 7.5	100 mM tri-Sodium citrate	7/B2
C3	20.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM Imidazole; pH 8.0	None	7/B3
C4	20.00 % w/v 2-Methyl-2,4-pentanediol	4.00 % w/v Glycerol	None	200 mM Sodium chloride	7/B4
C5	30.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM Sodium acetate; pH 4.6	20.00 mM Calcium chloride	7/B5
C6	30.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium acetate	7/B6
C7	30.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM HEPES; pH 7.5	500 mM Ammonium sulfate	7/C2
C8	30.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM HEPES; pH 7.5	200 mM tri-Sodium citrate	7/C3
C9	30.00 % w/v 2-Methyl-2,4-pentanediol	10.00 % w/v Polyethylene glycol 4,000	100 mM Imidazole; pH 8.0	None	7/C5
C10	30.00 % w/v 2-Methyl-2,4-pentanediol	20.00 % w/v Ethanol	None	None	7/C6
C11	35.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM Imidazole; pH 8.0	None	7/D2
C12	40.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM TRIS; pH 8.5	None	7/D3
D1	47.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM HEPES; pH 7.5	None	7/D4
D2	47.00 % w/v 2-Methyl-2,4-pentanediol	2.00 % w/v 2-Methyl-2-propanol	None	None	7/D5
D3	50.00 % w/v 2-Methyl-2,4-pentanediol	None	None	None	7/D6
D4	50.00 % w/v 2-Methyl-2,4-pentanediol	20.00 % w/v 2-Propanol	None	50.00 mM Sodium acetate, 50.00 mM Sodium chloride	8/A2
D5	50.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM TRIS; pH 8.5	100 mM Ammonium di-hydrogen phosphate	8/A3
D6	55.00 % w/v 2-Methyl-2,4-pentanediol	None	None	None	8/A4
D7	60.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM Sodium acetate; pH 4.6	10.00 mM Calcium chloride	8/A5
D8	70.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM MES; pH 6.5	None	8/B1
D9	70.00 % w/v 2-Methyl-2,4-pentanediol	None	100 mM TRIS; pH 8.5	None	8/B2
D10	20.00 % w/v Methanol	None	100 mM TRIS; pH 8.5	10.00 mM Calcium chloride	8/B3
D11	2.00 % w/v Ethanol	None	100 mM TRIS; pH 8.5	None	8/B4
D12	5.00 % w/v Ethanol	5.00 % w/v 2-Methyl-2,4-pentanediol	100 mM TRIS; pH 8.5	200 mM Sodium chloride	8/B6

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



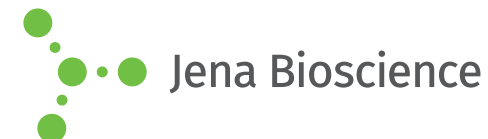


JBScreen Classic HTS II

(Ammonium Sulfate, MPD, Alcohol and Salt based)

Cat.-No.: CS-202L

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive	Classic Bulk No.
E1	10.00 % w/v Ethanol	None	100 mM TRIS; pH 8.5	None	8/C1
E2	12.00 % w/v Ethanol	4.00 % w/v Polyethylene glycol 400	100 mM Sodium acetate; pH 4.6	None	8/C2
E3	14.00 % w/v Ethanol	5.00 % w/v Glycerol	100 mM TRIS; pH 8.5	None	8/C3
E4	18.00 % w/v Ethanol	None	100 mM TRIS; pH 8.5	None	8/C4
E5	20.00 % w/v Ethanol	10.00 % w/v Glycerol	None	None	8/C6
E6	30.00 % w/v Ethanol	10.00 % w/v Polyethylene glycol 6,000	None	100 mM Sodium acetate	8/D1
E7	45.00 % w/v Ethanol	None	None	None	8/D2
E8	50.00 % w/v Ethanol	None	None	10.00 mM Sodium acetate	8/D3
E9	60.00 % w/v Ethanol	1.5 % w/v Polyethylene glycol 6,000	None	50.00 mM Sodium acetate	8/D4
E10	60.00 % w/v Ethanol	None	None	100 mM Sodium chloride	8/D5
E11	5.00 % w/v 2-Propanol	None	100 mM HEPES; pH 7.5	None	9/A1
E12	10.00 % w/v 2-Propanol	None	100 mM Sodium acetate; pH 4.6	200 mM Calcium chloride	9/A2
F1	10.00 % w/v 2-Propanol	None	100 mM HEPES; pH 7.5	200 mM tri-Sodium citrate	9/A3
F2	10.00 % w/v 2-Propanol	None	100 mM TRIS; pH 8.5	10.00 mM Magnesium chloride	9/A4
F3	15.00 % w/v 2-Propanol	None	100 mM MES; pH 6.5	200 mM tri-Sodium citrate	9/A6
F4	15.00 % w/v 2-Propanol	None	100 mM HEPES; pH 7.5	200 mM Magnesium chloride	9/B2
F5	15.00 % w/v 2-Propanol	None	100 mM TRIS; pH 8.5	200 mM Ammonium acetate	9/B3
F6	20.00 % w/v 2-Propanol	None	100 mM Sodium acetate; pH 4.6	200 mM Calcium chloride	9/B4
F7	25.00 % w/v 2-Propanol	None	100 mM HEPES; pH 7.5	100 mM Magnesium chloride	9/B6
F8	30.00 % w/v 2-Propanol	None	100 mM MES; pH 6.5	200 mM tri-Sodium citrate	9/C1
F9	30.00 % w/v 2-Propanol	None	100 mM TRIS; pH 8.5	200 mM Ammonium acetate	9/C3
F10	25.00 % w/v 2-Methyl-2-propanol	None	100 mM TRIS; pH 8.5	100 mM Calcium chloride	9/C4
F11	35.00 % w/v 2-Methyl-2-propanol	None	100 mM tri-Sodium citrate; pH 5.6	None	9/C5
F12	200 mM Ammonium di-hydrogen phosphate	None	None	None	9/C6

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





JBScreen Classic HTS II

(Ammonium Sulfate, MPD, Alcohol and Salt based)

Cat.-No.: CS-202L

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive	Classic Bulk No.
G1	200 mM Potassium Sodium tartrate	None	None	None	9/D1
G2	200 mM Magnesium acetate	None	None	None	9/D2
G3	400 mM Ammonium di-hydrogen phosphate	None	None	None	9/D3
G4	400 mM Potassium Sodium tartrate	None	100 mM TRIS; pH 8.5	None	9/D5
G5	500 mM Ammonium di-hydrogen phosphate	None	None	200 mM tri-Sodium citrate	9/D6
G6	500 mM Sodium acetate	None	100 mM Imidazole; pH 8.0	None	10/A1
G7	700 mM tri-Sodium citrate	None	100 mM HEPES; pH 7.5	None	10/A2
G8	800 mM Potassium Sodium tartrate	None	100 mM HEPES; pH 7.5	None	10/A4
G9	1.00 M Ammonium di-hydrogen phosphate	None	100 mM tri-Sodium citrate; pH 5.6	None	10/A5
G10	1.00 M Ammonium di-hydrogen phosphate	None	100 mM TRIS; pH 8.5	None	10/A6
G11	1.00 M Lithium sulfate	None	100 mM TRIS; pH 8.5	10.00 mM Nickel (II) chloride	10/B1
G12	1.00 M Sodium acetate	None	100 mM Imidazole; pH 8.0	None	10/B2
H1	1.4 M Sodium acetate	None	100 mM MES; pH 6.5	None	10/B4
H2	1.5 M Lithium sulfate	None	100 mM TRIS; pH 8.5	None	10/B6
H3	1.5 M tri-Sodium citrate; pH 6.5	None	None	None	10/C1
H4	1.6 M Magnesium sulfate	None	100 mM MES; pH 6.5	None	10/C2
H5	1.6 M Potassium Sodium tartrate	None	100 mM MES; pH 6.5	None	10/C3
H6	2.00 M Ammonium formate	None	100 mM MES; pH 6.5	None	10/C4
H7	2.00 M Ammonium di-hydrogen phosphate	None	100 mM TRIS; pH 8.5	None	10/C5
H8	2.00 M Sodium chloride	None	100 mM MES; pH 6.5	200 mM Sodium acetate	10/D2
H9	2.00 M Sodium formate	None	100 mM Sodium acetate; pH 4.6	None	10/D3
H10	1.00 M Ammonium di-hydrogen phosphate	30.00 % w/v Glycerol	100 mM TRIS; pH 8.5	None	10/D4
H11	4.00 M Sodium chloride	None	100 mM HEPES; pH 7.5	None	10/D5
H12	3.00 M Sodium formate	None	None	None	10/D6

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

