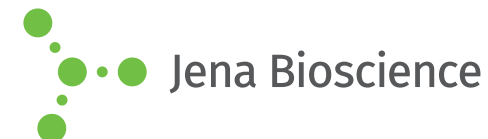




JBScreen Membrane HTS

Cat.-No.: CS-310

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
A1	15 % w/v Polyethylene glycol 400	15 % w/v Glycerol	100 mM HEPES; pH 7.5	200 mM Calcium chloride
A2	20 % w/v Polyethylene glycol 400	100 mM Sodium chloride	100 mM tri-Sodium citrate; pH 5.6	20 mM Magnesium chloride
A3	25 % w/v Polyethylene glycol 400	none	50 mM Sodium acetate; pH 4.6	50 mM Magnesium acetate
A4	30 % w/v Polyethylene glycol 400	50 mM Sodium sulfate	50 mM TRIS; pH 8.5	50 mM Lithium sulfate
A5	48 % w/v Polyethylene glycol 400	none	100 mM HEPES; pH 7.5	200 mM Calcium chloride
A6	20 % w/v Polyethylene glycol monomethyl ether 550	none	10 mM TRIS; pH 7.5	none
A7	30 % w/v Polyethylene glycol monomethyl ether 550	none	50 mM TRIS; pH 8.5	100 mM Magnesium chloride
A8	35 % w/v Polyethylene glycol 600	none	none	none
A9	28 % w/v Polyethylene glycol 1,000	10 % w/v Glycerol	100 mM TRICINE; pH 8.0	350 mM Sodium chloride
A10	10 % w/v Polyethylene glycol 1,500	5 % w/v Ethanol	none	100 mM Magnesium chloride, 100 mM Sodium chloride
A11	30 % w/v Polyethylene glycol 1,500	none	none	none
A12	5 % w/v Polyethylene glycol 2,000	none	none	none
B1	10 % w/v Polyethylene glycol 2,000	none	100 mM TRIS; pH 8.5	500 mM Magnesium chloride
B2	15 % w/v Polyethylene glycol 2,000	none	none	none
B3	15 % w/v Polyethylene glycol 2,000	none	none	100 mM Lithium chloride
B4	15 % w/v Polyethylene glycol 2,000	none	100 mM Sodium Phosphate; pH 6.2	20 mM tri-Sodium citrate
B5	15 % w/v Polyethylene glycol 2,000	none	100 mM Sodium Phosphate; pH 6.8	500 mM Sodium chloride
B6	15 % w/v Polyethylene glycol 2,000	none	20 mM BIS-TRIS; pH 7.0	none
B7	15 % w/v Polyethylene glycol 2,000	none	50 mM HEPES; pH 7.5	100 mM Magnesium chloride
B8	20 % w/v Polyethylene glycol 2,000	2 % w/v 2-Methyl-2,4-pentanediol	100 mM TRIS; pH 8.0	300 mM Magnesium nitrate
B9	25 % w/v Polyethylene glycol 2,000	15 % w/v Glycerol	100 mM BICINE; pH 9.0	300 mM Magnesium chloride
B10	30 % w/v Polyethylene glycol 2,000	none	200 mM Sodium Phosphate; pH 6.2	500 mM Sodium chloride
B11	8 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM Sodium acetate; pH 4.6	none
B12	10 % w/v Polyethylene glycol monomethyl ether 2,000	20 % w/v Glycerol	100 mM tri-Sodium citrate; pH 5.6	3 % w/v Polyethylene glycol 200

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





JBScreen Membrane HTS

Cat.-No.: CS-310

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
C1	12 % w/v Polyethylene glycol monomethyl ether 2,000	none	50 mM TRIS; pH 8.5	500 mM Sodium chloride
C2	10 % w/v Polyethylene glycol 3,350	none	50 mM tri-Sodium citrate; pH 5.6	150 mM Sodium chloride
C3	2 % w/v Polyethylene glycol 4,000	none	50 mM TRIS; pH 7.5	none
C4	5 % w/v Polyethylene glycol 4,000	none	none	none
C5	5 % w/v Polyethylene glycol 4,000	none	none	100 mM Potassium chloride
C6	5 % w/v Polyethylene glycol 4,000	10 % w/v Glycerol	50 mM MES; pH 6.5	100 mM Sodium chloride
C7	5 % w/v Polyethylene glycol 4,000	none	50 mM Sodium Phosphate; pH 6.7	none
C8	10 % w/v Polyethylene glycol 4,000	none	50 mM TRIS; pH 8.5	500 mM Sodium chloride
C9	12 % w/v Polyethylene glycol 4,000	none	100 mM ADA; pH 6.5	100 mM Lithium sulfate
C10	12 % w/v Polyethylene glycol 4,000	none	50 mM Sodium Phosphate; pH 6.8	none
C11	12 % w/v Polyethylene glycol 4,000	20 % w/v Glycerol	50 mM MOPS; pH 7.0	500 mM Potassium chloride
C12	15 % w/v Polyethylene glycol 4,000	none	10 mM TRIS; pH 7.5	100 mM Lithium chloride
D1	20 % w/v Polyethylene glycol 4,000	none	100 mM BIS-TRIS; pH 7.0	500 mM Sodium chloride
D2	20 % w/v Polyethylene glycol 4,000	none	100 mM Sodium Phosphate; pH 7.0	500 mM Sodium chloride
D3	20 % w/v Polyethylene glycol 4,000	150 mM Zinc acetate	50 mM TRIS; pH 7.5	50 mM Zinc chloride
D4	22 % w/v Polyethylene glycol 4,000	none	50 mM TRICINE; pH 8.0	none
D5	22 % w/v Polyethylene glycol 4,000	none	50 mM TRIS; pH 8.5	500 mM Sodium chloride
D6	30 % w/v Polyethylene glycol 4,000	none	none	none
D7	10 % w/v Polyethylene glycol monomethyl ether 5,000	none	100 mM tri-Sodium citrate; pH 5.6	100 mM Magnesium acetate
D8	5 % w/v Polyethylene glycol 6,000	none	none	100 mM Magnesium sulfate
D9	10 % w/v Polyethylene glycol 6,000	150 mM Zinc acetate	50 mM TRIS; pH 7.5	50 mM Zinc chloride
D10	15 % w/v Polyethylene glycol 6,000	none	50 mM di-Sodium succinate; pH 6.5	none
D11	12 % w/v Polyethylene glycol 8,000	10 % w/v 2-Methyl-2,4-pentanediol	none	25 mM Potassium di-hydrogen phosphate
D12	8 % w/v Polyethylene glycol 10,000	none	100 mM tri-Sodium citrate; pH 5.6	100 mM Magnesium acetate

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

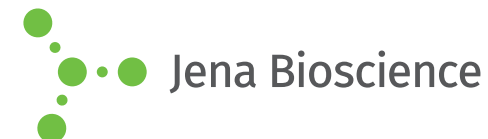




JBScreen Membrane HTS

Cat.-No.: CS-310

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
E1	500 mM Ammonium sulfate	none	1 M Sodium Potassium phosphate; pH 7.5	none
E2	1 M Ammonium sulfate	none	50 mM MES; pH 6.5	100 mM Zinc acetate
E3	1.2 M Ammonium sulfate	none	50 mM TRIS; pH 7.5	none
E4	1.2 M Ammonium sulfate	none	100 mM TRIS; pH 8.5	none
E5	1.4 M Ammonium sulfate	4 % w/v 2-Propanol	none	100 mM Ammonium acetate
E6	2 M Ammonium sulfate	none	none	none
E7	2 M Ammonium sulfate	none	100 mM tri-Sodium citrate; pH 5.6	none
E8	1 M Ammonium sulfate	2 % w/v Polyethylene glycol monomethyl ether 5,000	100 mM HEPES; pH 7.5	none
E9	3 M Ammonium sulfate	none	none	none
E10	3.5 M Ammonium sulfate	none	none	none
E11	3.5 M Ammonium sulfate	none	50 mM Sodium Potassium phosphate; pH 7.5	250 mM Sodium chloride
E12	25 % w/v 2-Methyl-2,4-pentanediol	none	100 mM BIS-TRIS; pH 7.0	none
F1	30 % w/v 2-Methyl-2,4-pentanediol	none	300 mM tri-Sodium citrate; pH 5.6	none
F2	25 % w/v Triethylene glycol	none	none	100 mM Glycine, 100 mM Ammonium sulfate
F3	30 % w/v 2-Propanol	20 % w/v Glycerol	100 mM Sodium acetate; pH 4.6	200 mM Calcium chloride
F4	none	none	50 mM Potassium Phosphate; pH 8.0	none
F5	none	none	100 mM tri-Sodium citrate; pH 4.8	none
F6	none	none	1 M Potassium Phosphate; pH 6.5	1 % w/v 1,4-Dioxane
F7	1 M tri-Sodium citrate	none	none	none
F8	1 M tri-Sodium citrate	none	none	500 mM Lithium chloride
F9	1.5 M Sodium chloride	none	100 mM Sodium acetate; pH 4.6	none
F10	none	none	1.5 M Potassium Phosphate; pH 7.0	none
F11	1.5 M Lithium sulfate	none	100 mM HEPES; pH 7.5	none
F12	2 M Sodium chloride	none	none	100 mM Sodium formate

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





JBScreen Membrane HTS

Cat.-No.: CS-310

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
G1	12 % w/v 2-Methyl-2,4-pentanediol	none	100 mM Sodium acetate; pH 4.6	100 mM Sodium chloride
G2	4 % w/v 2-Methyl-2,4-pentanediol	none	100 mM tri-Sodium citrate; pH 5.6	100 mM Magnesium chloride
G3	12 % w/v 2-Methyl-2,4-pentanediol	none	100 mM ADA; pH 6.5	none
G4	12 % w/v 2-Methyl-2,4-pentanediol	none	100 mM HEPES; pH 7.5	100 mM tri-Sodium citrate
G5	12 % w/v 2-Methyl-2,4-pentanediol	none	100 mM TRIS; pH 8.5	100 mM Lithium sulfate
G6	1 M tri-Sodium citrate	none	100 mM HEPES; pH 7.5	none
G7	200 mM Lithium sulfate	none	100 mM TRIS; pH 8.5	none
G8	100 mM Sodium chloride	none	100 mM tri-Sodium citrate; pH 5.6	none
G9	100 mM Sodium chloride	none	100 mM TRIS; pH 8.5	none
G10	1 M Ammonium di-hydrogen phosphate	none	100 mM Sodium acetate; pH 4.6	100 mM Lithium sulfate
G11	1 M Ammonium di-hydrogen phosphate	none	100 mM tri-Sodium citrate; pH 5.6	none
G12	1 M Ammonium di-hydrogen phosphate	none	100 mM ADA; pH 6.5	none
H1	2 M Ammonium di-hydrogen phosphate	none	100 mM TRIS; pH 7.5	none
H2	1 M Magnesium sulfate	none	100 mM Sodium acetate; pH 4.6	none
H3	1 M Magnesium sulfate	none	100 mM tri-Sodium citrate; pH 5.6	none
H4	1 M Magnesium sulfate	none	100 mM ADA; pH 6.5	100 mM Lithium sulfate
H5	400 mM Magnesium sulfate	none	50 mM TRIS; pH 7.5	none
H6	400 mM Magnesium sulfate	none	100 mM TRIS; pH 8.5	100 mM Potassium Sodium tartrate
H7	1 M Potassium Sodium tartrate	none	100 mM HEPES; pH 7.5	none
H8	100 mM Potassium Sodium tartrate	none	100 mM HEPES; pH 7.5	100 mM Lithium sulfate
H9	4 M Sodium formate	none	none	none
H10	2 M Sodium formate; pH 4.6	none	none	none
H11	1.4 M Sodium acetate	none	100 mM MES; pH 6.5	none
H12	100 mM Sodium acetate	none	100 mM TRIS; pH 8.5	none

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

