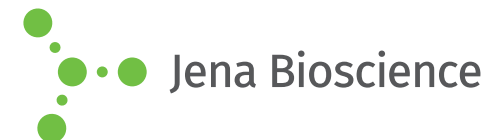




JBScreen Nuc-Pro HTS

Cat.-No.: CS-209L

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
A1	20 % v/v Polyethylene glycol 200	none	50 mM HEPES; pH 7.5	200 mM Potassium chloride, 25 mM Magnesium sulfate
A2	50 % v/v Polyethylene glycol 200	none	100 mM TRIS; pH 8.0	none
A3	5 % v/v Polyethylene glycol 400	none	50 mM PIPES; pH 7.0	30 mM Magnesium chloride
A4	5 % v/v Polyethylene glycol 400	none	20 mM MES; pH 5.8	15 mM Magnesium formate, 2 mM Cobalt (II) chloride
A5	10 % v/v Polyethylene glycol 400	none	50 mM HEPES; pH 7.0	100 mM Potassium chloride
A6	15 % v/v Polyethylene glycol 400	none	50 mM MES; pH 6.5	80 mM Magnesium acetate, 15 mM Magnesium chloride
A7	15 % v/v Polyethylene glycol 400	none	100 mM TRIS; pH 8.0	80 mM Calcium chloride, 20 mM Sodium chloride
A8	20 % v/v Polyethylene glycol 400	none	50 mM BIS-TRIS propane; pH 6.8	60 mM Magnesium chloride
A9	25 % v/v Polyethylene glycol 400	none	100 mM tri-Sodium citrate; pH 5.6	130 mM Sodium chloride, 60 mM Magnesium chloride
A10	30 % v/v Polyethylene glycol 400	none	100 mM HEPES; pH 7.5	200 mM Calcium chloride
A11	30 % v/v Polyethylene glycol 400	none	50 mM TRIS; pH 8.5	100 mM Potassium chloride, 10 mM Magnesium chloride
A12	25 % v/v Polyethylene glycol monomethyl ether 550	none	50 mM HEPES; pH 7.0	10 mM Magnesium chloride
B1	20 % w/v Polyethylene glycol 1,000	none	50 mM MES; pH 6.5	200 mM Magnesium chloride, 100 mM Sodium chloride
B2	30 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate
B3	10 % w/v Polyethylene glycol 3,350	none	100 mM MES; pH 6.5	100 mM Calcium chloride, 13 % v/v Glycerol
B4	25 % w/v Polyethylene glycol 3,350	none	50 mM MES; pH 6.0	200 mM Sodium formate, 10 % v/v Glycerol
B5	30 % w/v Polyethylene glycol 3,350	none	50 mM di-Sodium succinate; pH 5.5	100 mM Sodium chloride
B6	35 % w/v Polyethylene glycol 3,350	none	50 mM TRIS; pH 7.5	50 mM Potassium chloride
B7	5 % w/v Polyethylene glycol 4,000	none	50 mM HEPES; pH 7.0	200 mM Ammonium sulfate, 20 mM Magnesium acetate
B8	5 % w/v Polyethylene glycol 4,000	none	50 mM MES; pH 6.0	5 mM Magnesium sulfate
B9	10 % w/v Polyethylene glycol 4,000	none	50 mM MES; pH 6.5	200 mM Ammonium acetate, 10 mM Calcium chloride
B10	10 % w/v Polyethylene glycol 4,000	none	50 mM Imidazole; pH 7.2	20 mM Zinc sulfate
B11	15 % w/v Polyethylene glycol 4,000	none	50 mM TRIS; pH 7.5	150 mM Potassium chloride, 20 mM Magnesium chloride
B12	15 % w/v Polyethylene glycol 4,000	none	50 mM tri-Sodium citrate; pH 5.0	100 mM Sodium chloride, 20 mM Ammonium sulfate

*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
C1	20 % w/v Polyethylene glycol 4,000	none	100 mM tri-Sodium citrate; pH 5.6	100 mM Potassium chloride, 10 % v/v 2-Propanol
C2	20 % w/v Polyethylene glycol 4,000	none	50 mM MOPS; pH 7.0	100 mM Sodium chloride
C3	20 % w/v Polyethylene glycol 4,000	none	100 mM HEPES; pH 7.5	none
C4	24 % w/v Polyethylene glycol 4,000	none	50 mM di-Sodium succinate; pH 5.5	60 mM Magnesium chloride, 300 mM Sodium chloride
C5	25 % w/v Polyethylene glycol 4,000	none	50 mM tri-Sodium citrate; pH 5.0	50 mM Ammonium sulfate
C6	30 % w/v Polyethylene glycol 4,000	none	50 mM MES; pH 6.5	80 mM Magnesium acetate
C7	30 % w/v Polyethylene glycol 4,000	none	50 mM TRIS; pH 8.5	150 mM Ammonium chloride, 10 mM Calcium chloride
C8	32 % w/v Polyethylene glycol 4,000	none	100 mM TRIS; pH 8.5	5 % v/v Glycerol
C9	36 % w/v Polyethylene glycol 4,000	none	50 mM Sodium acetate; pH 5.0	none
C10	5 % w/v Polyethylene glycol 6,000	none	20 mM BIS-TRIS; pH 6.0	60 mM di-Ammonium hydrogen citrate, 20 mM Magnesium chloride
C11	10 % w/v Polyethylene glycol 6,000	none	50 mM tri-Sodium citrate; pH 4.0	100 mM Sodium chloride
C12	10 % w/v Polyethylene glycol 6,000	none	50 mM HEPES; pH 7.0	200 mM Ammonium acetate, 150 mM Magnesium acetate
D1	15 % w/v Polyethylene glycol 6,000	none	10 mM TRIS; pH 7.5	none
D2	18 % w/v Polyethylene glycol 6,000	none	50 mM Sodium acetate; pH 5.0	none
D3	20 % w/v Polyethylene glycol 6,000	none	50 mM BIS-TRIS propane; pH 7.0	7 % v/v 2-Methyl-2,4-pentanediol, 5 % v/v 2-Methyl-2-propanol
D4	20 % w/v Polyethylene glycol 6,000	none	50 mM di-Potassium L-malate; pH 5.0	30 mM Calcium chloride
D5	5 % w/v Polyethylene glycol 8,000	none	50 mM HEPES; pH 7.5	20 mM Magnesium chloride
D6	10 % w/v Polyethylene glycol 8,000	none	100 mM TRIS; pH 8.0	10 % v/v Glycerol, 1 mM Tris(2-carboxyethyl)phosphine hydrochloride
D7	10 % w/v Polyethylene glycol 8,000	none	50 mM MES; pH 6.5	200 mM Potassium chloride, 100 mM Magnesium acetate
D8	15 % w/v Polyethylene glycol 8,000	none	50 mM BIS-TRIS propane; pH 6.8	100 mM Ammonium sulfate, 10 % v/v Glycerol
D9	15 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Calcium acetate
D10	18 % w/v Polyethylene glycol 8,000	none	100 mM TRIS; pH 8.0	200 mM Magnesium formate
D11	20 % w/v Polyethylene glycol 8,000	none	10 mM TRIS; pH 7.5	10 mM Calcium chloride
D12	15 % w/v Polyethylene glycol 20,000	none	100 mM MES; pH 6.5	60 mM Manganese (II) 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
E1	1.2 M Ammonium sulfate	none	100 mM HEPES; pH 7.5	2 % w/v Polyethylene glycol 400
E2	20 mM Calcium chloride	none	100 mM Glycine; pH 8.0	none
E3	2 M Lithium chloride	none	50 mM MES; pH 6.0	200 mM Calcium acetate, 1 mM Cobalt (II) chloride
E4	600 mM Lithium sulfate	none	50 mM MES; pH 6.0	10 mM Magnesium chloride
E5	1 M Lithium sulfate	8 % w/v Polyethylene glycol 400	50 mM HEPES; pH 7.5	none
E6	1 M Lithium sulfate	none	none	50 mM tri-Sodium citrate, 3 % w/v 2-Propanol
E7	1.2 M Lithium sulfate	none	50 mM MES; pH 6.5	30 mM Magnesium chloride
E8	1.2 M Lithium sulfate	none	50 mM MES; pH 6.5	50 mM Magnesium chloride, 2 mM Cobalt (II) chloride
E9	1.5 M Lithium sulfate	none	50 mM TRIS; pH 8.5	5 % w/v Glycerol
E10	1.6 M Lithium sulfate	2 % w/v Polyethylene glycol 1,000	50 mM HEPES; pH 7.5	none
E11	1.7 M Lithium sulfate	none	50 mM HEPES; pH 7.0	50 mM Magnesium sulfate
E12	1.7 M Lithium sulfate	none	none	10 % w/v Glycerol
F1	2 M Lithium sulfate	none	none	3 % w/v 2-Methyl-2,4-pentanediol
F2	5 mM Magnesium chloride	none	50 mM MES; pH 6.5	2.5 mM Cobalt (II) chloride
F3	40 mM Magnesium chloride	none	50 mM MES; pH 6.0	none
F4	80 mM Magnesium chloride	none	50 mM HEPES; pH 7.5	none
F5	200 mM Sodium chloride	none	50 mM BIS-TRIS propane; pH 7.0	none
F6	400 mM Sodium chloride	none	50 mM BIS-TRIS propane; pH 6.8	none
F7	600 mM Sodium chloride	none	50 mM MES; pH 6.0	100 mM Ammonium acetate, 5 mM Magnesium sulfate
F8	2.5 M Sodium chloride	none	50 mM TRIS; pH 7.5	200 mM Magnesium chloride
F9	1.8 M Sodium formate	none	100 mM TRIS; pH 8.0	none
F10	2 M Sodium formate	none	100 mM Sodium acetate; pH 4.6	none
F11	100 mM di-Sodium hydrogen phosphate	none	none	80 mM Sodium chloride
F12	1 M di-Sodium tartrate	none	50 mM TRIS; pH 7.5	30 mM Magnesium chloride

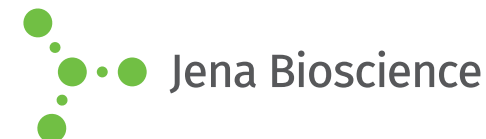
*pH values indicated are those of the 1.0 M buffer stock solution prior to dilution with other components



JBScreen Nuc-Pro HTS

Cat.-No.: CS-209L

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
G1	10 % w/v 1,6-Hexanediol	none	50 mM MES; pH 6.5	20 mM Magnesium chloride
G2	20 % w/v 1,6-Hexanediol	none	50 mM HEPES; pH 7.0	50 mM Ammonium chloride, 10 mM Magnesium chloride
G3	35 % w/v 1,6-Hexanediol	none	50 mM TRIS; pH 8.5	75 mM Magnesium sulfate
G4	10 % v/v 2-Methyl-2,4-pentanediol	none	50 mM TRIS; pH 7.5	50 mM Ammonium acetate
G5	10 % v/v 2-Methyl-2,4-pentanediol	none	50 mM HEPES; pH 7.0	80 mM Potassium chloride, 10 mM Magnesium sulfate
G6	15 % v/v 2-Methyl-2,4-pentanediol	none	50 mM ADA; pH 6.5	100 mM Sodium acetate
G7	15 % v/v 2-Methyl-2,4-pentanediol	none	50 mM di-Sodium succinate; pH 5.5	10 mM Magnesium acetate
G8	18 % v/v 2-Methyl-2,4-pentanediol	none	20 mM MES; pH 5.8	10 mM Magnesium chloride
G9	23 % v/v 2-Methyl-2,4-pentanediol	none	50 mM MES; pH 6.0	100 mM Sodium chloride
G10	26 % v/v 2-Methyl-2,4-pentanediol	none	50 mM PIPES; pH 7.0	65 mM Magnesium chloride, 1 mM Cobalt (III) Hexamine chloride
G11	27 % v/v 2-Methyl-2,4-pentanediol	none	20 mM MES; pH 5.8	400 mM Sodium chloride, 120 mM Calcium chloride
G12	35 % v/v 2-Methyl-2,4-pentanediol	none	20 mM BIS-TRIS; pH 6.0	50 mM Sodium chloride, 10 mM Calcium chloride
H1	50 % v/v 2-Methyl-2,4-pentanediol	none	100 mM tri-Sodium citrate; pH 5.6	10 mM Magnesium chloride
H2	5 % v/v 2-Propanol	none	50 mM TRIS; pH 7.5	10 mM Magnesium chloride
H3	5 % v/v 2-Propanol	none	50 mM MES; pH 6.5	100 mM Calcium acetate
H4	9 % v/v 2-Propanol	none	50 mM Imidazole; pH 7.2	15 mM Magnesium acetate, 15 mM Magnesium chloride
H5	10 % v/v 2-Propanol	none	50 mM MES; pH 6.5	80 mM Ammonium acetate
H6	10 % v/v 2-Propanol	none	50 mM di-Sodium succinate; pH 5.5	2 mM Cobalt (II) chloride
H7	13 % v/v 2-Propanol	none	50 mM MOPS; pH 7.0	200 mM Potassium chloride, 6 mM Cobalt (III) Hexamine chloride
H8	15 % v/v 2-Propanol	none	50 mM MES; pH 6.0	20 mM Magnesium chloride
H9	10 % v/v 1,4-Dioxane	none	50 mM HEPES; pH 7.5	none
H10	10 % v/v Ethanol	none	50 mM MES; pH 6.5	20 mM Magnesium chloride, 1 mM Cobalt (II) chloride
H11	20 % v/v Ethylene glycol	5 % w/v Polyethylene glycol 3,350	none	20 mM Magnesium chloride
H12	15 % v/v Glycerol	none	100 mM Sodium acetate; pH 4.6	200 mM Sodium chloride

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

