



Cat.-No.: CS-213L

SCREEN FORMULATION



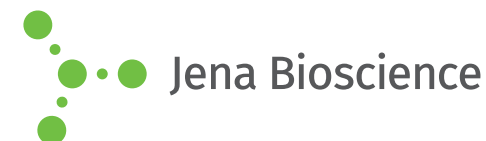
No.	Precipitant	Buffer	Additive
A1	15 % v/v Polyethylene glycol 300	100 mM HEPES; pH 7.6	50 mM Ethylenediaminetetraacetic acid disodium salt; pH 8.0, 1.2 % v/v 1,7 Heptanediol
A2	30 % v/v Polyethylene glycol 300	100 mM HEPES; pH 7.5	100 mM Ammonium di-hydrogen phosphate, 2 % v/v 2-Methyl-2,4-pentanediol
A3	30 % v/v Polyethylene glycol 300	100 mM Sodium Phosphate; pH 6.3	150 mM Sodium chloride
A4	32 % v/v Polyethylene glycol 300	100 mM HEPES; pH 7.5	100 mM Ammonium di-hydrogen phosphate, 1 % v/v 1,7 Heptanediol
A5	37 % v/v Polyethylene glycol 300	100 mM BIS-TRIS propane; pH 6.5	100 mM di-Ammonium hydrogen phosphate
A6	10 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	150 mM Ammonium sulfate
A7	13 % v/v Polyethylene glycol 400	100 mM MES; pH 6.0	170 mM Potassium Sodium tartrate, 0.45 % v/v Jeffamine® M-600; pH 7.0
A8	14 % v/v Polyethylene glycol 400	100 mM TRIS; pH 7.5	6 % v/v 2-Methyl-2,4-pentanediol
A9	18 % v/v Polyethylene glycol 400	100 mM TRIS; pH 7.5	100 mM tri-Sodium citrate
A10	18 % v/v Polyethylene glycol 400	50 mM HEPES; pH 7.0	100 mM Potassium sulfate
A11	18.5 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 5.0	50 mM Lithium sulfate, 100 mM Sodium chloride
A12	19.5 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.0	350 mM Ammonium di-hydrogen phosphate
B1	20 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 5.5	300 mM di-Sodium malonate, 5 mM 2-Aminoethanesulfonic acid
B2	20 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.0	400 mM Sodium chloride
B3	20 % v/v Polyethylene glycol 400	100 mM MES; pH 6.5	400 mM Potassium nitrate; pH 6.9, 1 mM Tris(2-carboxyethyl)phosphine hydrochloride
B4	21 % v/v Polyethylene glycol 400	100 mM MES; pH 6.5	70 mM di-Ammonium hydrogen phosphate
B5	23 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 5.25	300 mM di-Sodium malonate, 5 mM Nickel (II) chloride
B6	25 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	200 mM Potassium Sodium tartrate
B7	26 % v/v Polyethylene glycol 400	100 mM MES; pH 6.7	200 mM Lithium sulfate, 3.5 % v/v 1,4-Butanediol, 4 % v/v Dimethyl sulfoxide
B8	26 % v/v Polyethylene glycol 400	100 mM MES; pH 6.0	300 mM di-Sodium malonate, 5 mM Strontium chloride
B9	26 % v/v Polyethylene glycol 400	100 mM TRIS; pH 8.0	300 mM Ammonium sulfate
B10	26.5 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 5.0	50 mM Sodium thiocyanate, 2 % v/v 2,5-Hexanediol
B11	27 % v/v Polyethylene glycol 400	100 mM TRIS; pH 7.75	220 mM Sodium formate, 5 % v/v 1,4-Butanediol
B12	27 % v/v Polyethylene glycol 400	100 mM HEPES; pH 6.9	150 mM Ammonium fluoride, 2.5 % v/v Jeffamine® M-600

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



Cat.-No.: CS-213L

SCREEN FORMULATION



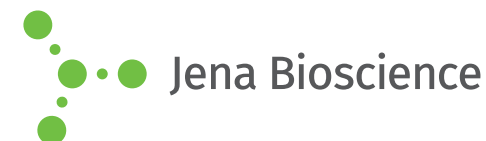
No.	Precipitant	Buffer	Additive
C1	27.5 % v/v Polyethylene glycol 400	100 mM BIS-TRIS propane; pH 6.7	120 mM di-Sodium tartrate, 3 % v/v 1,3-Butanediol
C2	27.5 % v/v Polyethylene glycol 400	100 mM BIS-TRIS propane; pH 6.4	150 mM Potassium Sodium tartrate
C3	28 % v/v Polyethylene glycol 400	100 mM BIS-TRIS propane; pH 7.0	300 mM Potassium formate
C4	28 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 4.5	300 mM Ammonium di-hydrogen phosphate, 10 mM Magnesium chloride
C5	28.5 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.0	100 mM di-Ammonium hydrogen phosphate, 6 mM Tris(2-carboxyethyl)phosphine hydrochloride
C6	29 % v/v Polyethylene glycol 400	100 mM MES; pH 6.8	200 mM di-Ammonium hydrogen phosphate
C7	29 % v/v Polyethylene glycol 400	50 mM tri-Sodium citrate; pH 4.0	200 mM Lithium sulfate
C8	29.5 % v/v Polyethylene glycol 400	100 mM TRIS; pH 7.75	350 mM Sodium formate, 5 % v/v 1,4-Butanediol
C9	30 % v/v Polyethylene glycol 400	100 mM TRIS; pH 8.0	200 mM di-Sodium malonate
C10	30 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 6.5	185 mM Lithium sulfate
C11	30 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 5.0	200 mM Magnesium chloride
C12	30 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 5.5	200 mM Lithium sulfate
D1	30 % v/v Polyethylene glycol 400	100 mM BIS-TRIS propane; pH 7.5	300 mM Ammonium acetate, 2 % w/v D-(+)-Glucose
D2	30 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.0	100 mM Sodium chloride
D3	30 % v/v Polyethylene glycol 400	100 mM MES; pH 6.0	100 mM Magnesium sulfate, 2.5 % v/v Polypropylene glycol 400
D4	30 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 6.0	400 mM Potassium nitrate
D5	30 % v/v Polyethylene glycol 400	100 mM TRIS; pH 7.5	400 mM Lithium chloride
D6	30 % v/v Polyethylene glycol 400	100 mM TRIS; pH 8.0	100 mM Magnesium sulfate
D7	30 % v/v Polyethylene glycol 400	100 mM MES; pH 6.3	100 mM Ammonium formate
D8	30 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.0	100 mM Sodium thiocyanate
D9	30 % v/v Polyethylene glycol 400	100 mM MES; pH 6.5	100 mM Sodium thiocyanate, 20 mM Calcium chloride
D10	30 % v/v Polyethylene glycol 400	100 mM TRIS; pH 8.0	100 mM Sodium thiocyanate, 20 mM Calcium chloride
D11	30 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 5.5	100 mM Sodium chloride, 3 % w/v D-(+)-Trehalose
D12	30 % v/v Polyethylene glycol 400	50 mM MES; pH 6.5	100 mM Magnesium chloride

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



Cat.-No.: CS-213L

SCREEN FORMULATION



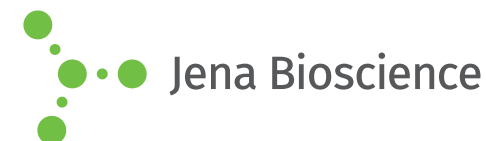
No.	Precipitant	Buffer	Additive
E1	30 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	2 mM β -Mercaptoethanol
E2	31 % v/v Polyethylene glycol 400	100 mM MES; pH 6.0	300 mM di-Sodium malonate, 5 mM Cobalt (III) Hexamine chloride
E3	31 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	150 mM tri-Sodium citrate, 350 mM Magnesium chloride
E4	32 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.8	70 mM Ammonium fluoride, 6 % v/v Polypropylene glycol 400
E5	32 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.0	100 mM Sodium chloride
E6	32 % v/v Polyethylene glycol 400	50 mM MES; pH 6.5	100 mM Sodium thiocyanate, 20 mM Calcium chloride
E7	32 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 4.5	75 mM Sodium chloride, 130 mM Magnesium chloride
E8	32.5 % v/v Polyethylene glycol 400	100 mM BIS-TRIS propane; pH 6.75	150 mM Sodium sulfate, 6 % v/v 1,4-Butanediol
E9	32.5 % v/v Polyethylene glycol 400	100 mM MES; pH 6.2	100 mM Potassium Sodium tartrate, 5 % v/v Ethylene glycol
E10	34 % v/v Polyethylene glycol 400	100 mM TRIS; pH 8.7	90 mM tri-Sodium citrate, 120 mM Ammonium sulfate
E11	34 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.0	300 mM Lithium sulfate, 7.5 % v/v Dimethyl sulfoxide
E12	35 % v/v Polyethylene glycol 400	none	4 % v/v Glycerol, 80 mM tri-Sodium citrate, 100 mM TRICINE
F1	35 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 5.0	370 mM Ammonium acetate, 3 % v/v 1-Propanol
F2	36 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.2	100 mM Ammonium chloride
F3	37.5 % v/v Polyethylene glycol 400	100 mM tri-Sodium citrate; pH 6.0	180 mM di-Ammonium tartrate, 4 % v/v 2-Methyl-2,4-pentanediol
F4	38 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.0	300 mM Lithium chloride, 30 mM Strontium chloride
F5	39 % v/v Polyethylene glycol 400	100 mM TRIS; pH 8.5	100 mM Potassium chloride
F6	39.8 % v/v Polyethylene glycol 400	100 mM BIS-TRIS propane; pH 7.2	100 mM Ammonium di-hydrogen phosphate
F7	40 % v/v Polyethylene glycol 400	100 mM TRIS; pH 8.0	4 % v/v Dimethyl sulfoxide, 1 % v/v 1,7 Heptanediol
F8	40 % v/v Polyethylene glycol 400	50 mM MES; pH 6.5	1.6 M Sodium chloride
F9	40 % v/v Polyethylene glycol 400	100 mM ADA; pH 6.5	160 mM Lithium chloride, 4 mM Strontium chloride
F10	40 % v/v Polyethylene glycol 400	100 mM ADA; pH 6.5	200 mM Lithium sulfate
F11	42 % v/v Polyethylene glycol 400	100 mM MES; pH 6.5	150 mM Sodium acetate
F12	25 % v/v Polyethylene glycol dimethyl ether 500	100 mM MES; pH 6.0	10 mM Copper (II) chloride, 200 mM Ammonium formate

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



Cat.-No.: CS-213L

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
G1	30 % v/v Polyethylene glycol dimethyl ether 500	100 mM MES; pH 6.0	2 mM Cadmium chloride
G2	30 % v/v Polyethylene glycol dimethyl ether 500	100 mM tri-Sodium citrate; pH 6.0	100 mM Magnesium chloride, 100 mM Sodium chloride, 100 mM Ammonium sulfate
G3	14 % v/v Polyethylene glycol monomethyl ether 550	100 mM HEPES; pH 7.5	200 mM Lithium sulfate
G4	15 % v/v Polyethylene glycol monomethyl ether 550	100 mM HEPES; pH 7.0	200 mM Potassium Sodium tartrate, 1 mM Tris(2-carboxyethyl)phosphine hydrochloride
G5	25 % v/v Polyethylene glycol monomethyl ether 550	50 mM ADA; pH 6.25	350 mM Sodium nitrate, 50 mM di-Sodium malonate; pH 7.0
G6	28 % v/v Polyethylene glycol monomethyl ether 550	100 mM TRIS; pH 8.0	100 mM Lithium sulfate
G7	28 % v/v Polyethylene glycol monomethyl ether 550	50 mM ADA; pH 7.0	250 mM Ammonium sulfate
G8	25 % v/v Polyethylene glycol 600	100 mM ADA; pH 7.0	none
G9	22.5 % w/v Polyethylene glycol 1,500	100 mM MES; pH 5.5	300 mM Sodium acetate
G10	12 % w/v Polyethylene glycol 3,350	100 mM Sodium acetate; pH 4.6	200 mM di-Sodium malonate
G11	25 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS; pH 5.5	200 mM Lithium sulfate
G12	25 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS; pH 5.5	none
H1	10 % w/v Polyethylene glycol 4,000	100 mM Sodium acetate; pH 4.6	200 mM Potassium chloride
H2	20 % w/v Polyethylene glycol 6,000	100 mM BICINE; pH 9.0	none
H3	30 % w/v Polyethylene glycol 8,000	none	200 mM Ammonium sulfate
H4	10 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM MES; pH 6.5	100 mM Ammonium chloride, 10 mM Calcium chloride
H5	20 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM TRIS; pH 8.0	100 mM Potassium formate
H6	4 % v/v 2-Methyl-2,4-pentanediol	100 mM tri-Sodium citrate; pH 5.6	100 mM Sodium chloride, 100 mM Lithium nitrate
H7	5 % v/v 2-Methyl-2,4-pentanediol	100 mM tri-Sodium citrate; pH 5.6	100 mM Sodium chloride, 60 mM Magnesium acetate
H8	8 % v/v 2-Methyl-2,4-pentanediol	100 mM ADA; pH 6.7	400 mM Potassium nitrate, 100 mM tri-Potassium citrate
H9	none	1 M Sodium Potassium phosphate; pH 5.1	300 mM D-(+)-Trehalose
H10	1.5 M Sodium chloride	75 mM Sodium acetate; pH 4.6	none
H11	1 M Sodium acetate	100 mM MES; pH 6.5	none
H12	1 M Lithium sulfate	100 mM TRIS; pH 8.5	10 mM Nickel (II) chloride

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