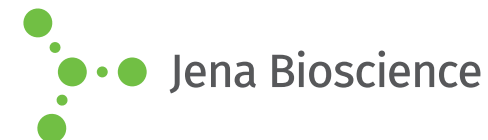




Cat.-No.: CS-128

SCREEN FORMULATION



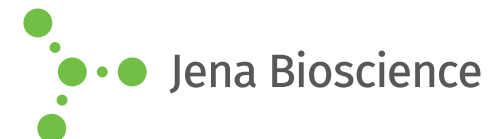
No.	Precipitant	Buffer	Additive
A1	22.5 % v/v Polyethylene glycol 200	50 mM Sodium acetate; pH 4.8	none
A2	20 % v/v Polyethylene glycol 300	50 mM Sodium acetate; pH 5.2	none
A3	20 % v/v Polyethylene glycol monomethyl ether 350	50 mM MES; pH 5.6	none
A4	20 % v/v Polyethylene glycol 400	50 mM MES; pH 6.0	none
A5	20 % v/v Polyethylene glycol monomethyl ether 550	50 mM ADA; pH 6.4	none
A6	20 % v/v Polyethylene glycol 600	50 mM ADA; pH 6.8	none
A7	17.5 % w/v Polyethylene glycol 1,000	50 mM HEPES; pH 7.1	none
A8	17.5 % w/v Polyethylene glycol 1,500	50 mM HEPES; pH 7.3	none
A9	15 % w/v Polyethylene glycol 2,000	50 mM TRIS; pH 7.6	none
A10	15 % w/v Polyethylene glycol monomethyl ether 2,000	50 mM TRIS; pH 8.0	none
A11	15 % w/v Polyethylene glycol 3,000	50 mM BICINE; pH 8.4	none
A12	15 % w/v Polyethylene glycol 4,000	50 mM BICINE; pH 8.8	none
B1	15 % w/v Polyethylene glycol 1,500	50 mM Sodium acetate; pH 4.8	3.6 % w/v Polyethylene glycol 4,000
B2	12.9 % w/v Polyethylene glycol 2,000	50 mM Sodium acetate; pH 5.2	6.4 % v/v Polyethylene glycol 200
B3	12.9 % w/v Polyethylene glycol monomethyl ether 2,000	50 mM MES; pH 5.6	5.7 % v/v Polyethylene glycol 300
B4	12.9 % w/v Polyethylene glycol 3,000	50 mM MES; pH 6.0	5.7 % v/v Polyethylene glycol monomethyl ether 350
B5	12.9 % w/v Polyethylene glycol 4,000	50 mM ADA; pH 6.4	5.7 % v/v Polyethylene glycol 400
B6	19.3 % v/v Polyethylene glycol 200	50 mM ADA; pH 6.8	5.7 % v/v Polyethylene glycol monomethyl ether 550
B7	17.1 % v/v Polyethylene glycol 300	50 mM HEPES; pH 7.1	5.7 % v/v Polyethylene glycol 600
B8	17.1 % v/v Polyethylene glycol monomethyl ether 350	50 mM HEPES; pH 7.3	5 % w/v Polyethylene glycol 1,000
B9	17.1 % v/v Polyethylene glycol 400	50 mM TRIS; pH 7.6	5 % w/v Polyethylene glycol 1,500
B10	17.1 % v/v Polyethylene glycol monomethyl ether 550	50 mM TRIS; pH 8.0	4.3 % w/v Polyethylene glycol 2,000
B11	17.1 % v/v Polyethylene glycol 600	50 mM BICINE; pH 8.4	4.3 % w/v Polyethylene glycol monomethyl ether 2,000
B12	15 % w/v Polyethylene glycol 1,000	50 mM BICINE; pH 8.8	4.3 % w/v Polyethylene glycol 3,000

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



Cat.-No.: CS-128

SCREEN FORMULATION



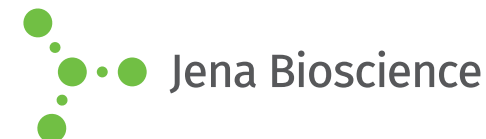
No.	Precipitant	Buffer	Additive
C1	14.3 % v/v Polyethylene glycol monomethyl ether 350	50 mM Sodium acetate; pH 4.8	11.4 % v/v Polyethylene glycol 300
C2	14.3 % v/v Polyethylene glycol 400	50 mM Sodium acetate; pH 5.2	11.4 % v/v Polyethylene glycol monomethyl ether 350
C3	14.3 % v/v Polyethylene glycol monomethyl ether 550	50 mM MES; pH 5.6	11.4 % v/v Polyethylene glycol 400
C4	14.3 % v/v Polyethylene glycol 600	50 mM MES; pH 6.0	11.4 % v/v Polyethylene glycol monomethyl ether 550
C5	12.5 % w/v Polyethylene glycol 1,000	50 mM ADA; pH 6.4	11.4 % v/v Polyethylene glycol 600
C6	12.5 % w/v Polyethylene glycol 1,500	50 mM ADA; pH 6.8	10 % w/v Polyethylene glycol 1,000
C7	10.7 % w/v Polyethylene glycol 2,000	50 mM HEPES; pH 7.1	10 % w/v Polyethylene glycol 1,500
C8	10.7 % w/v Polyethylene glycol monomethyl ether 2,000	50 mM HEPES; pH 7.3	8.6 % w/v Polyethylene glycol 2,000
C9	10.7 % w/v Polyethylene glycol 3,000	50 mM TRIS; pH 7.6	8.6 % w/v Polyethylene glycol monomethyl ether 2,000
C10	10.7 % w/v Polyethylene glycol 4,000	50 mM TRIS; pH 8.0	8.6 % w/v Polyethylene glycol 3,000
C11	16.1 % v/v Polyethylene glycol 200	50 mM BICINE; pH 8.4	7.1 % w/v Polyethylene glycol 4,000
C12	14.3 % v/v Polyethylene glycol 300	50 mM BICINE; pH 8.8	12.9 % v/v Polyethylene glycol 200
D1	8.6 % w/v Polyethylene glycol monomethyl ether 2,000	50 mM Sodium acetate; pH 4.8	17.1 % v/v Polyethylene glycol 400
D2	8.6 % w/v Polyethylene glycol 3,000	50 mM Sodium acetate; pH 5.2	17.1 % v/v Polyethylene glycol monomethyl ether 550
D3	8.6 % w/v Polyethylene glycol 4,000	50 mM MES; pH 5.6	17.1 % v/v Polyethylene glycol 600
D4	12.9 % v/v Polyethylene glycol 200	50 mM MES; pH 6.0	15 % w/v Polyethylene glycol 1,000
D5	11.4 % v/v Polyethylene glycol 300	50 mM ADA; pH 6.4	15 % w/v Polyethylene glycol 1,500
D6	11.4 % v/v Polyethylene glycol monomethyl ether 350	50 mM ADA; pH 6.8	12.9 % w/v Polyethylene glycol 2,000
D7	11.4 % v/v Polyethylene glycol 400	50 mM HEPES; pH 7.1	12.9 % w/v Polyethylene glycol monomethyl ether 2,000
D8	11.4 % v/v Polyethylene glycol monomethyl ether 550	50 mM HEPES; pH 7.3	12.9 % w/v Polyethylene glycol 3,000
D9	11.4 % v/v Polyethylene glycol 600	50 mM TRIS; pH 7.6	10.7 % w/v Polyethylene glycol 4,000
D10	10 % w/v Polyethylene glycol 1,000	50 mM TRIS; pH 8.0	19.3 % v/v Polyethylene glycol 200
D11	10 % w/v Polyethylene glycol 1,500	50 mM BICINE; pH 8.4	17.1 % v/v Polyethylene glycol 300
D12	8.6 % w/v Polyethylene glycol 2,000	50 mM BICINE; pH 8.8	17.1 % v/v Polyethylene glycol monomethyl ether 350

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



Cat.-No.: CS-128

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
E1	8.6 % v/v Polyethylene glycol monomethyl ether 550	50 mM Sodium acetate; pH 4.8	22.9 % v/v Polyethylene glycol 600
E2	8.6 % v/v Polyethylene glycol 600	50 mM Sodium acetate; pH 5.2	20 % w/v Polyethylene glycol 1,000
E3	7.5 % w/v Polyethylene glycol 1,000	50 mM MES; pH 5.6	20 % w/v Polyethylene glycol 1,500
E4	7.5 % w/v Polyethylene glycol 1,500	50 mM MES; pH 6.0	17.1 % w/v Polyethylene glycol 2,000
E5	6.4 % w/v Polyethylene glycol 2,000	50 mM ADA; pH 6.4	17.1 % w/v Polyethylene glycol monomethyl ether 2,000
E6	6.4 % w/v Polyethylene glycol monomethyl ether 2,000	50 mM ADA; pH 6.8	17.1 % w/v Polyethylene glycol 3,000
E7	6.4 % w/v Polyethylene glycol 3,000	50 mM HEPES; pH 7.1	14.3 % w/v Polyethylene glycol 4,000
E8	6.4 % w/v Polyethylene glycol 4,000	50 mM HEPES; pH 7.3	25.7 % v/v Polyethylene glycol 200
E9	9.6 % v/v Polyethylene glycol 200	50 mM TRIS; pH 7.6	22.9 % v/v Polyethylene glycol 300
E10	8.6 % v/v Polyethylene glycol 300	50 mM TRIS; pH 8.0	22.9 % v/v Polyethylene glycol monomethyl ether 350
E11	8.6 % v/v Polyethylene glycol monomethyl ether 350	50 mM BICINE; pH 8.4	22.9 % v/v Polyethylene glycol 400
E12	8.6 % v/v Polyethylene glycol 400	50 mM BICINE; pH 8.8	22.9 % v/v Polyethylene glycol monomethyl ether 550
F1	4.3 % w/v Polyethylene glycol 4,000	50 mM Sodium acetate; pH 4.8	25 % w/v Polyethylene glycol 1,500
F2	6.4 % v/v Polyethylene glycol 200	50 mM Sodium acetate; pH 5.2	21.4 % w/v Polyethylene glycol 2,000
F3	5.7 % v/v Polyethylene glycol 300	50 mM MES; pH 5.6	21.4 % w/v Polyethylene glycol monomethyl ether 2,000
F4	5.7 % v/v Polyethylene glycol monomethyl ether 350	50 mM MES; pH 6.0	21.4 % w/v Polyethylene glycol 3,000
F5	5.7 % v/v Polyethylene glycol 400	50 mM ADA; pH 6.4	17.9 % w/v Polyethylene glycol 4,000
F6	5.7 % v/v Polyethylene glycol monomethyl ether 550	50 mM ADA; pH 6.8	32.1 % v/v Polyethylene glycol 200
F7	5.7 % v/v Polyethylene glycol 600	50 mM HEPES; pH 7.1	28.6 % v/v Polyethylene glycol 300
F8	5 % w/v Polyethylene glycol 1,000	50 mM HEPES; pH 7.3	28.6 % v/v Polyethylene glycol monomethyl ether 350
F9	5 % w/v Polyethylene glycol 1,500	50 mM TRIS; pH 7.6	28.6 % v/v Polyethylene glycol 400
F10	4.3 % w/v Polyethylene glycol 2,000	50 mM TRIS; pH 8.0	28.6 % v/v Polyethylene glycol monomethyl ether 550
F11	4.3 % w/v Polyethylene glycol monomethyl ether 2,000	50 mM BICINE; pH 8.4	28.6 % v/v Polyethylene glycol 600
F12	4.3 % w/v Polyethylene glycol 3,000	50 mM BICINE; pH 8.8	25 % w/v Polyethylene glycol 1,000

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



Cat.-No.: CS-128

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
G1	2.5 % w/v Polyethylene glycol 1,000	50 mM Sodium acetate; pH 4.8	25.7 % w/v Polyethylene glycol monomethyl ether 2,000
G2	2.5 % w/v Polyethylene glycol 1,500	50 mM Sodium acetate; pH 5.2	25.7 % w/v Polyethylene glycol 3,000
G3	2.1 % w/v Polyethylene glycol 2,000	50 mM MES; pH 5.6	21.4 % w/v Polyethylene glycol 4,000
G4	2.1 % w/v Polyethylene glycol monomethyl ether 2,000	50 mM MES; pH 6.0	38.6 % v/v Polyethylene glycol 200
G5	2.1 % w/v Polyethylene glycol 3,000	50 mM ADA; pH 6.4	34.3 % v/v Polyethylene glycol 300
G6	2.1 % w/v Polyethylene glycol 4,000	50 mM ADA; pH 6.8	34.3 % v/v Polyethylene glycol monomethyl ether 350
G7	3.2 % v/v Polyethylene glycol 200	50 mM HEPES; pH 7.1	34.3 % v/v Polyethylene glycol 400
G8	2.9 % v/v Polyethylene glycol 300	50 mM HEPES; pH 7.3	34.3 % v/v Polyethylene glycol monomethyl ether 550
G9	2.9 % v/v Polyethylene glycol monomethyl ether 350	50 mM TRIS; pH 7.6	34.3 % v/v Polyethylene glycol 600
G10	2.9 % v/v Polyethylene glycol 400	50 mM TRIS; pH 8.0	30 % w/v Polyethylene glycol 1,000
G11	2.9 % v/v Polyethylene glycol monomethyl ether 550	50 mM BICINE; pH 8.4	30 % w/v Polyethylene glycol 1,500
G12	2.9 % v/v Polyethylene glycol 600	50 mM BICINE; pH 8.8	25.7 % w/v Polyethylene glycol 2,000
H1	none	50 mM Sodium acetate; pH 4.8	25 % w/v Polyethylene glycol 4,000
H2	none	50 mM Sodium acetate; pH 5.2	45 % v/v Polyethylene glycol 200
H3	none	50 mM MES; pH 5.6	40 % v/v Polyethylene glycol 300
H4	none	50 mM MES; pH 6.0	40 % v/v Polyethylene glycol monomethyl ether 350
H5	none	50 mM ADA; pH 6.4	40 % v/v Polyethylene glycol 400
H6	none	50 mM ADA; pH 6.8	40 % v/v Polyethylene glycol monomethyl ether 550
H7	none	50 mM HEPES; pH 7.1	40 % v/v Polyethylene glycol 600
H8	none	50 mM HEPES; pH 7.3	35 % w/v Polyethylene glycol 1,000
H9	none	50 mM TRIS; pH 7.6	35 % w/v Polyethylene glycol 1,500
H10	none	50 mM TRIS; pH 8.0	30 % w/v Polyethylene glycol 2,000
H11	none	50 mM BICINE; pH 8.4	30 % w/v Polyethylene glycol monomethyl ether 2,000
H12	none	50 mM BICINE; pH 8.8	30 % w/v Polyethylene glycol 3,000

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