



JBScreen Kinase HTS

Cat.-No.: CS-204L

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
A1	1 M Ammonium sulfate	100 mM tri-Sodium citrate; pH 5.6	200 mM Magnesium acetate, 10 mM Dithiothreitol
A2	1.3 M Ammonium sulfate	100 mM tri-Sodium citrate; pH 5.6	none
A3	1.3 M Ammonium sulfate	100 mM TRIS; pH 8.5	none
A4	1.8 M Ammonium sulfate	100 mM MES; pH 6.5	25 mM Cobalt (II) chloride
A5	2 M Ammonium sulfate	100 mM tri-Sodium citrate; pH 3.1	200 mM Sodium chloride
A6	2 M Ammonium sulfate	100 mM Sodium acetate; pH 4.6	50 mM Magnesium chloride
A7	2 M Ammonium sulfate	100 mM Sodium acetate; pH 4.6	none
A8	2 M Ammonium sulfate	100 mM HEPES; pH 7.5	2 % v/v Polyethylene glycol monomethyl ether 550
A9	2 M Ammonium sulfate	100 mM HEPES; pH 7.5	none
A10	2 M Ammonium sulfate	100 mM TRIS; pH 8.5	6 mM Magnesium chloride
A11	1.5 M Lithium sulfate	100 mM TRIS; pH 8.5	10 mM Nickel sulfate
A12	1 M Lithium chloride	100 mM tri-Sodium citrate; pH 4.2	none
B1	2 M Sodium chloride	100 mM Sodium acetate; pH 4.6	none
B2	2 M Sodium chloride	100 mM MES; pH 6.5	100 mM di-Sodium hydrogen phosphate, 100 mM Potassium di-hydrogen phosphate
B3	3.3 M Sodium chloride	100 mM HEPES; pH 7.5	1 % v/v Glycerol
B4	1.2 M Sodium acetate	100 mM MES; pH 6.5	6.3 mM Calcium chloride
B5	3.7 M Sodium formate	100 mM BICINE; pH 9.5	2 % w/v Polyethylene glycol 3,000
B6	500 mM di-Sodium malonate; pH 6.0	50 mM PIPES; pH 6.0	1.6 % v/v Glycerol, 10 mM Dithiothreitol
B7	500 mM di-Sodium hydrogen phosphate	100 mM CAPS; pH 10.0	500 mM Potassium di-hydrogen phosphate, 200 mM Lithium sulfate
B8	1.2 M di-Sodium tartrate	100 mM TRIS; pH 8.5	5 mM Dithiothreitol
B9	1 M Potassium Sodium tartrate	100 mM MES; pH 6.5	none
B10	30 % v/v Jeffamine® M-600	100 mM MES; pH 6.5	50 mM Cesium chloride
B11	40 % v/v 2-Methyl-2,4-pentanediol	100 mM MES; pH 6.5	none
B12	50 % v/v 2-Methyl-2,4-pentanediol	100 mM HEPES; pH 7.5	none

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





JBScreen Kinase HTS

Cat.-No.: CS-204L

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
C1	10 % v/v Polyethylene glycol 400	50 mM TRIS; pH 8.5	1 mM Dithiothreitol, 1 mM Ethylenediaminetetraacetic acid disodium salt, 300 mM Sodium chloride
C2	15 % v/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	200 mM Calcium chloride
C3	25 % v/v Polyethylene glycol 400	100 mM MES; pH 6.5	10 % v/v 2-Propanol
C4	25 % v/v Polyethylene glycol 400	100 mM TRIS; pH 8.5	150 mM tri-Sodium citrate
C5	15 % v/v Polyethylene glycol monomethyl ether 550	100 mM Sodium acetate; pH 4.6	5 % v/v Ethylene glycol
C6	20 % v/v Polyethylene glycol monomethyl ether 550	100 mM BICINE; pH 9.0	100 mM Sodium chloride
C7	20 % w/v Polyethylene glycol 1,000	100 mM TRIS; pH 8.5	1 mM Dithiothreitol
C8	35 % w/v Polyethylene glycol 1,000	100 mM HEPES; pH 7.5	50 mM Lithium sulfate
C9	12 % w/v Polyethylene glycol 2,000	100 mM MES; pH 6.5	200 mM Magnesium acetate
C10	25 % w/v Polyethylene glycol 2,000	100 mM Sodium acetate; pH 4.6	100 mM Magnesium chloride
C11	30 % w/v Polyethylene glycol 2,000	100 mM Sodium acetate; pH 4.6	50 mM Magnesium chloride
C12	24 % w/v Polyethylene glycol monomethyl ether 2,000	100 mM Potassium phosphate citrate; pH 5.0	none
D1	12 % w/v Polyethylene glycol 3,350	100 mM MES; pH 6.5	500 mM Sodium chloride
D2	12 % w/v Polyethylene glycol 3,350	50 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium sulfate, 50 mM Magnesium sulfate
D3	15 % w/v Polyethylene glycol 3,350	100 mM Imidazole; pH 7.5	250 mM Ammonium sulfate, 10 mM Cadmium chloride
D4	20 % w/v Polyethylene glycol 3,350	150 mM di-Sodium DL-malate; pH 7.0	none
D5	20 % w/v Polyethylene glycol 3,350	100 mM HEPES; pH 7.5	200 mM Sodium chloride, 20 mM L-Glutathione reduced
D6	20 % w/v Polyethylene glycol 3,350	100 mM TRIS; pH 8.5	120 mM Sodium chloride, 5 mM Dithiothreitol
D7	20 % w/v Polyethylene glycol 3,350	none	200 mM Potassium nitrate
D8	22 % w/v Polyethylene glycol 3,350	none	100 mM Ammonium formate
D9	24 % w/v Polyethylene glycol 3,350	100 mM tri-Sodium citrate; pH 5.0	none
D10	30 % w/v Polyethylene glycol 3,350	100 mM Sodium acetate; pH 4.6	200 mM Ammonium acetate
D11	30 % w/v Polyethylene glycol 3,350	200 mM Ammonium acetate; pH 5.6	20 % v/v 2-Propanol, 200 mM Calcium chloride
D12	32.5 % w/v Polyethylene glycol 3,350	100 mM TRIS; pH 8.5	500 mM Sodium chloride, 200 mM Magnesium chloride

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

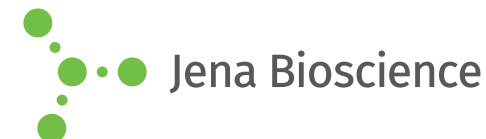




JBScreen Kinase HTS

Cat.-No.: CS-204L

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
E1	8 % w/v Polyethylene glycol 4,000	50 mM MES; pH 6.5	10 mM Magnesium chloride, 10 mM Dithiothreitol
E2	10 % w/v Polyethylene glycol 4,000	50 mM PIPES; pH 7.0	10 mM Dithiothreitol
E3	10 % w/v Polyethylene glycol 4,000	100 mM HEPES; pH 7.5	15 % v/v Ethylene glycol, 10 % v/v 2-Propanol, 200 mM Magnesium chloride
E4	10 % w/v Polyethylene glycol 4,000	100 mM HEPES; pH 7.5	none
E5	15 % w/v Polyethylene glycol 4,000	100 mM HEPES; pH 7.5	10 % v/v 2-Propanol
E6	15 % w/v Polyethylene glycol 4,000	75 mM TRIS; pH 8.5	200 mM Sodium chloride, 1 % w/v Polyethylene glycol 6,000, 75 mM Sodium acetate
E7	15 % w/v Polyethylene glycol 4,000	100 mM di-Sodium DL-malate; pH 5.5	200 mM Ammonium sulfate
E8	20 % w/v Polyethylene glycol 4,000	100 mM BIS-TRIS; pH 6.5	100 mM Sodium chloride
E9	20 % w/v Polyethylene glycol 4,000	100 mM HEPES; pH 7.5	200 mM Magnesium chloride
E10	20 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Magnesium chloride
E11	25 % w/v Polyethylene glycol 4,000	100 mM MES; pH 6.5	200 mM Magnesium chloride
E12	25 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	100 mM Lithium sulfate
F1	28 % w/v Polyethylene glycol 4,000	200 mM Lithium acetate; pH 7.5	none
F2	30 % w/v Polyethylene glycol 4,000	100 mM MES; pH 6.5	200 mM Sodium acetate
F3	30 % w/v Polyethylene glycol 4,000	150 mM TRIS; pH 8.5	200 mM Ammonium sulfate
F4	8 % w/v Polyethylene glycol monomethyl ether 5,000	100 mM HEPES; pH 7.5	10 % v/v 2-Propanol
F5	25 % w/v Polyethylene glycol monomethyl ether 5,000	100 mM MES; pH 6.5	200 mM Ammonium sulfate
F6	30 % w/v Polyethylene glycol monomethyl ether 5,000	100 mM HEPES; pH 7.5	200 mM Ammonium sulfate
F7	30 % w/v Polyethylene glycol monomethyl ether 5,000	100 mM ADA; pH 6.5	100 mM Ammonium sulfate
F8	20 % w/v Polyethylene glycol 6,000	100 mM MES; pH 6.5	none
F9	28 % w/v Polyethylene glycol 6,000	100 mM MES; pH 6.5	10 mM Dithiothreitol
F10	30 % w/v Polyethylene glycol 6,000	100 mM HEPES; pH 7.5	175 mM Lithium sulfate
F11	30 % w/v Polyethylene glycol 6,000	100 mM PIPES; pH 7.0	10 mM Dithiothreitol
F12	32 % w/v Polyethylene glycol 6,000	100 mM MES; pH 6.5	none

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

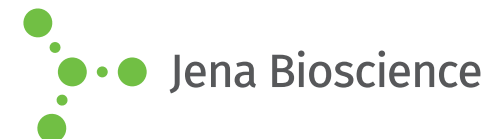




JBScreen Kinase HTS

Cat.-No.: CS-204L

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
G1	7 % w/v Polyethylene glycol 8,000	100 mM MES; pH 6.5	20 % v/v Ethylene glycol
G2	7 % w/v Polyethylene glycol 8,000	100 mM MES; pH 6.5	150 mM Calcium acetate, 16 % v/v Ethylene glycol
G3	10 % w/v Polyethylene glycol 8,000	100 mM TRIS; pH 8.5	10 % v/v Polyethylene glycol 200
G4	12 % w/v Polyethylene glycol 8,000	100 mM HEPES; pH 7.5	none
G5	12 % w/v Polyethylene glycol 8,000	100 mM TRIS; pH 8.5	250 mM di-Sodium tartrate
G6	16 % w/v Polyethylene glycol 8,000	100 mM HEPES; pH 7.5	100 mM Potassium di-hydrogen phosphate
G7	16 % w/v Polyethylene glycol 8,000	100 mM HEPES; pH 7.5	150 mM Sodium chloride, 2 % v/v Ethylene glycol
G8	18 % w/v Polyethylene glycol 8,000	100 mM MES; pH 6.5	200 mM Magnesium acetate
G9	18 % w/v Polyethylene glycol 8,000	100 mM MES; pH 6.5	none
G10	18 % w/v Polyethylene glycol 8,000	100 mM TRIS; pH 8.5	none
G11	20 % w/v Polyethylene glycol 8,000	100 mM tri-Sodium citrate; pH 5.0	100 mM Magnesium acetate
G12	20 % w/v Polyethylene glycol 8,000	100 mM TRIS; pH 8.5	200 mM Magnesium chloride, 2 % v/v Ethylene glycol
H1	22 % w/v Polyethylene glycol 8,000	100 mM TRIS; pH 8.5	2 % v/v Ethylene glycol
H2	25 % w/v Polyethylene glycol 8,000	100 mM Sodium acetate; pH 4.6	50 mM Magnesium chloride
H3	30 % w/v Polyethylene glycol 8,000	100 mM MES; pH 6.5	200 mM Ammonium sulfate, 4 % v/v 1,3-Propanediol
H4	30 % w/v Polyethylene glycol 8,000	100 mM HEPES; pH 7.5	10 mM Dithiothreitol, 20 % v/v Glycerol
H5	9 % w/v Polyethylene glycol 8,000	100 mM MES; pH 6.5	200 mM Zinc acetate
H6	16 % w/v Polyethylene glycol 10,000	100 mM BIS-TRIS; pH 6.5	300 mM Ammonium sulfate, 5 % v/v Ethylene glycol
H7	10 % w/v Polyethylene glycol 10,000	100 mM HEPES; pH 7.5	8 % v/v Ethylene glycol
H8	15 % w/v Polyethylene glycol 10,000	100 mM HEPES; pH 7.5	5 mM Dithiothreitol
H9	15 % w/v Polyethylene glycol 10,000	100 mM TRIS; pH 8.5	none
H10	12 % w/v Polyethylene glycol 20,000	100 mM MES; pH 6.5	none
H11	10 % w/v Polyethylene glycol 20,000	100 mM HEPES; pH 7.5	100 mM Ammonium formate
H12	15 % w/v Polyethylene glycol 20,000	none	10 mM Potassium hydrogen tartrate

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

