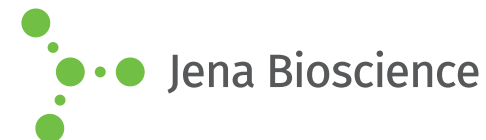




JBScreen Kinase 1

Cat.-No.: CS-131

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	6.25 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

