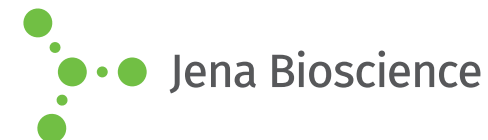




Cat.-No.: CS-164

## SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
G1	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM Sodium fluoride
G2	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM Sodium bromide
G3	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM Sodium iodide
G4	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM Potassium thiocyanate
G5	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM Sodium nitrate
G6	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM Sodium formate
G7	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM Sodium acetate
G8	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM Sodium sulfate
G9	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM Potassium Sodium tartrate
G10	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	none
G11	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM tri-Sodium citrate
G12	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 7.5	200 mM di-Sodium malonate
H1	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM Sodium fluoride
H2	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM Sodium bromide
H3	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM Sodium iodide
H4	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM Potassium thiocyanate
H5	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM Sodium nitrate
H6	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM Sodium formate
H7	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM Sodium acetate
H8	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM Sodium sulfate
H9	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM Potassium Sodium tartrate
H10	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	20 mM Sodium Potassium phosphate
H11	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM tri-Sodium citrate
H12	20 % w/v Polyethylene glycol 3,350	100 mM BIS-TRIS propane; pH 8.5	200 mM di-Sodium malonate

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