



JBScreen Pentaerythritol 3

(PEE 270 based)

Cat.-No.: CS-193

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
E1	25 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM Sodium acetate; pH 4.6	none
E2	35 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM Sodium acetate; pH 4.6	none
E3	45 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM Sodium acetate; pH 4.6	none
E4	25 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM MES; pH 6.5	none
E5	35 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM MES; pH 6.5	none
E6	45 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM MES; pH 6.5	none
E7	25 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM HEPES; pH 7.5	none
E8	35 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM HEPES; pH 7.5	none
E9	45 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM HEPES; pH 7.5	none
E10	25 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM TRIS; pH 8.5	none
E11	35 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM TRIS; pH 8.5	none
E12	45 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM TRIS; pH 8.5	none
F1	25 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM Sodium acetate; pH 4.6	50 mM Magnesium chloride
F2	35 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate
F3	45 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM Sodium acetate; pH 4.6	400 mM Potassium chloride
F4	25 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM MES; pH 6.5	50 mM Magnesium chloride
F5	35 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM MES; pH 6.5	200 mM Ammonium sulfate
F6	45 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM MES; pH 6.5	400 mM Potassium chloride
F7	25 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM HEPES; pH 7.5	50 mM Magnesium chloride
F8	35 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM HEPES; pH 7.5	200 mM Ammonium sulfate
F9	45 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM HEPES; pH 7.5	400 mM Potassium chloride
F10	25 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM TRIS; pH 8.5	50 mM Magnesium chloride
F11	35 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM TRIS; pH 8.5	200 mM Ammonium sulfate
F12	45 % w/v Pentaerythritol ethoxylate (3/4 EO/OH)	100 mM TRIS; pH 8.5	400 mM Potassium chloride

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

