



## JBScreen Pentaerythritol 2

(PEP 629 based)

Cat.-No.: CS-192

## SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
C1	25 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM Sodium acetate; pH 4.6	none
C2	35 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM Sodium acetate; pH 4.6	none
C3	45 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM Sodium acetate; pH 4.6	none
C4	25 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM MES; pH 6.5	none
C5	35 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM MES; pH 6.5	none
C6	45 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM MES; pH 6.5	none
C7	25 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM HEPES; pH 7.5	none
C8	35 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM HEPES; pH 7.5	none
C9	45 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM HEPES; pH 7.5	none
C10	25 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM TRIS; pH 8.5	none
C11	35 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM TRIS; pH 8.5	none
C12	45 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM TRIS; pH 8.5	none
D1	25 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM Sodium acetate; pH 4.6	50 mM Magnesium chloride
D2	35 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate
D3	45 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM Sodium acetate; pH 4.6	300 mM Potassium chloride
D4	25 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM MES; pH 6.5	50 mM Magnesium chloride
D5	35 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM MES; pH 6.5	200 mM Ammonium sulfate
D6	45 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM MES; pH 6.5	300 mM Potassium chloride
D7	25 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM HEPES; pH 7.5	50 mM Magnesium chloride
D8	35 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM HEPES; pH 7.5	200 mM Ammonium sulfate
D9	45 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM HEPES; pH 7.5	300 mM Potassium chloride
D10	25 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM TRIS; pH 8.5	50 mM Magnesium chloride
D11	35 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM TRIS; pH 8.5	200 mM Ammonium sulfate
D12	45 % w/v Pentaerythritol propoxylate (17/8 PO/OH)	100 mM TRIS; pH 8.5	300 mM Potassium chloride

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

