



JBScreen Pentaerythritol HTS

Cat.-No.: CS-210L

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
A1	25 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM Sodium acetate; pH 4.6	none
A2	35 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM Sodium acetate; pH 4.6	none
A3	45 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM Sodium acetate; pH 4.6	none
A4	25 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM MES; pH 6.5	none
A5	35 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM MES; pH 6.5	none
A6	45 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM MES; pH 6.5	none
A7	25 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM HEPES; pH 7.5	none
A8	35 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM HEPES; pH 7.5	none
A9	45 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM HEPES; pH 7.5	none
A10	25 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM TRIS; pH 8.5	none
A11	35 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM TRIS; pH 8.5	none
A12	45 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM TRIS; pH 8.5	none
B1	25 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM Sodium acetate; pH 4.6	50 mM Magnesium chloride
B2	35 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate
B3	45 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM Sodium acetate; pH 4.6	400 mM Potassium chloride
B4	25 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM MES; pH 6.5	50 mM Magnesium chloride
B5	35 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM MES; pH 6.5	200 mM Ammonium sulfate
B6	45 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM MES; pH 6.5	400 mM Potassium chloride
B7	25 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM HEPES; pH 7.5	50 mM Magnesium chloride
B8	35 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM HEPES; pH 7.5	200 mM Ammonium sulfate
B9	45 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM HEPES; pH 7.5	400 mM Potassium chloride
B10	25 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM TRIS; pH 8.5	50 mM Magnesium chloride
B11	35 % w/v Pentaerythritol Propoxylate (5/4 PO/OH)	100 mM TRIS; pH 8.5	200 mM Ammonium sulfate
B12	45 % w/v Pentaerythritol Propoxylate (5/4 PO/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

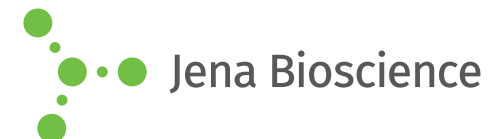




JBScreen Pentaerythritol HTS

Cat.-No.: CS-210L

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

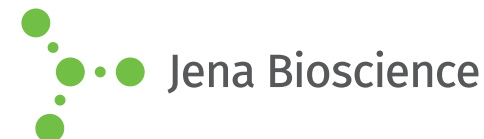




JBScreen Pentaerythritol HTS

Cat.-No.: CS-210L

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





JBScreen Pentaerythritol HTS

Cat.-No.: CS-210L

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
G1	25 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM Sodium acetate; pH 4.6	none
G2	35 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM Sodium acetate; pH 4.6	none
G3	45 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM Sodium acetate; pH 4.6	none
G4	25 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM MES; pH 6.5	none
G5	35 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM MES; pH 6.5	none
G6	45 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM MES; pH 6.5	none
G7	25 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM HEPES; pH 7.5	none
G8	35 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM HEPES; pH 7.5	none
G9	45 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM HEPES; pH 7.5	none
G10	25 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM TRIS; pH 8.5	none
G11	35 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM TRIS; pH 8.5	none
G12	45 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM TRIS; pH 8.5	none
H1	25 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM Sodium acetate; pH 4.6	50 mM Magnesium chloride
H2	35 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate
H3	45 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM Sodium acetate; pH 4.6	400 mM Potassium chloride
H4	25 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM MES; pH 6.5	50 mM Magnesium chloride
H5	35 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM MES; pH 6.5	200 mM Ammonium sulfate
H6	45 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM MES; pH 6.5	400 mM Potassium chloride
H7	25 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM HEPES; pH 7.5	50 mM Magnesium chloride
H8	35 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM HEPES; pH 7.5	200 mM Ammonium sulfate
H9	45 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM HEPES; pH 7.5	400 mM Potassium chloride
H10	25 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM TRIS; pH 8.5	50 mM Magnesium chloride
H11	35 % w/v Pentaerythritol ethoxylate (15/4 EO/OH)	100 mM TRIS; pH 8.5	200 mM Ammonium sulfate
H12	45 % w/v Pentaerythritol ethoxylate (15/4 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

