

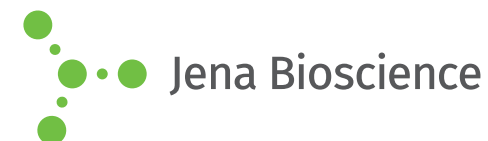


JBScreen Pentaerythritol 2

(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

