



JBScreen Classic 8

(MPD/Alcohol based)

Cat.-No.: CS-108L

SCREEN FORMULATION



Jena Bioscience

| No. | Precipitant 1 | Precipitant 2 | Buffer | Additive |
|------|-----------------------------------|-------------------------------------|-------------------------------|---|
| 8/A1 | 50 % w/v 2-Methyl-2,4-pentanediol | 15 % w/v Ethanol | none | 10 mM Sodium acetate |
| 8/A2 | 50 % w/v 2-Methyl-2,4-pentanediol | 20 % w/v 2-Propanol | none | 50 mM Sodium acetate, 50 mM Sodium chloride |
| 8/A3 | 50 % w/v 2-Methyl-2,4-pentanediol | none | 100 mM TRIS; pH 8.5 | 100 mM Ammonium di-hydrogen phosphate |
| 8/A4 | 55 % w/v 2-Methyl-2,4-pentanediol | none | none | none |
| 8/A5 | 60 % w/v 2-Methyl-2,4-pentanediol | none | 100 mM Sodium acetate; pH 4.6 | 10 mM Calcium chloride |
| 8/A6 | 60 % w/v 2-Methyl-2,4-pentanediol | none | none | 20 mM Sodium acetate |
| 8/B1 | 70 % w/v 2-Methyl-2,4-pentanediol | none | 100 mM MES; pH 6.5 | none |
| 8/B2 | 70 % w/v 2-Methyl-2,4-pentanediol | none | 100 mM TRIS; pH 8.5 | none |
| 8/B3 | 20 % w/v Methanol | none | 100 mM TRIS; pH 8.5 | 10 mM Calcium chloride |
| 8/B4 | 2 % w/v Ethanol | none | 100 mM TRIS; pH 8.5 | none |
| 8/B5 | 5 % w/v Ethanol | 5 % w/v 2-Methyl-2,4-pentanediol | 100 mM HEPES; pH 7.5 | none |
| 8/B6 | 5 % w/v Ethanol | 5 % w/v 2-Methyl-2,4-pentanediol | 100 mM TRIS; pH 8.5 | 200 mM Sodium chloride |
| 8/C1 | 10 % w/v Ethanol | none | 100 mM TRIS; pH 8.5 | none |
| 8/C2 | 12 % w/v Ethanol | 4 % w/v Polyethylene glycol 400 | 100 mM Sodium acetate; pH 4.6 | none |
| 8/C3 | 14 % w/v Ethanol | 5 % w/v Glycerol | 100 mM TRIS; pH 8.5 | none |
| 8/C4 | 18 % w/v Ethanol | none | 100 mM TRIS; pH 8.5 | none |
| 8/C5 | 20 % w/v Ethanol | none | none | none |
| 8/C6 | 20 % w/v Ethanol | 10 % w/v Glycerol | none | none |
| 8/D1 | 30 % w/v Ethanol | 10 % w/v Polyethylene glycol 6,000 | none | 100 mM Sodium acetate |
| 8/D2 | 45 % w/v Ethanol | none | none | none |
| 8/D3 | 50 % w/v Ethanol | none | none | 10 mM Sodium acetate |
| 8/D4 | 60 % w/v Ethanol | 1.5 % w/v Polyethylene glycol 6,000 | none | 50 mM Sodium acetate |
| 8/D5 | 60 % w/v Ethanol | none | none | 100 mM Sodium chloride |
| 8/D6 | 2 % w/v 2-Propanol | none | 100 mM TRIS; pH 8.5 | 10 mM Magnesium sulfate |

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

