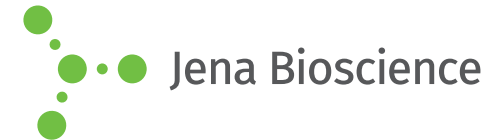




JBScreen Plus Chaotropic

Cat.-No.: CS-502

SCREEN FORMULATION



Jena Bioscience

| No. | Additive | Type | CAS Number |
|-----|-------------------------------------|-------------------|------------|
| A1 | 1.00 M Urea | Non ionic | 57-13-6 |
| A2 | 4.00 M Dimethyl sulfoxide | Non ionic | 67-68-5 |
| A3 | 4.00 M Dimethylformamide | Non ionic | 68-12-2 |
| A4 | 1.00 M Pyridine | Non ionic | 110-86-1 |
| A5 | 0.50 M e-Caprolactam | Non ionic | 105-60-2 |
| A6 | 0.10 M Phenol | Non ionic | 108-95-2 |
| B1 | 2.00 M Trimethylamine hydrochloride | Chaotropic Cation | 593-81-7 |
| B2 | 2.00 M Guanidine hydrochloride | Chaotropic Cation | 50-01-1 |
| B3 | 2.00 M Tetramethylammonium Chloride | Chaotropic Cation | 75-57-0 |
| B4 | 1.00 M Cesium chloride | Chaotropic Cation | 7647-17-8 |
| B5 | 1.00 M Rubidium chloride | Chaotropic Cation | 7791-11-9 |
| B6 | 1.00 M Potassium chloride | Chaotropic Cation | 7447-40-7 |
| C1 | 2.00 M Potassium thiocyanate | Chaotropic Anion | 333-20-0 |
| C2 | 2.00 M Potassium Cyanate | Chaotropic Anion | 590-28-3 |
| C3 | 1.00 M Potassium nitrate | Chaotropic Anion | 7757-79-1 |
| C4 | 1.00 M Potassium acetate | Chaotropic Anion | 127-08-2 |
| C5 | 1.00 M Potassium iodide | Chaotropic Anion | 7681-11-0 |
| C6 | 1.00 M Potassium bromide | Chaotropic Anion | 7758-02-3 |
| D1 | 2.00 M Sodium trichloroacetate | Chaotropic Anion | 650-51-1 |
| D2 | 2.00 M Sodium thiocyanate | Chaotropic Anion | 540-72-7 |
| D3 | 2.00 M Sodium perchlorate | Chaotropic Anion | 7601-89-0 |
| D4 | 2.00 M Sodium nitrate | Chaotropic Anion | 7631-99-4 |
| D5 | 1.00 M Sodium iodide | Chaotropic Anion | 7681-82-5 |
| D6 | 2.00 M Sodium bromide | Chaotropic Anion | 7647-15-6 |

