

JBScreen Cryo 1

Cat.-No: CC-103

| Number | Cryoprotectant 1 | Cryoprotectant 2 | Cryoprotectant 3 | Buffer | pH | Additive 1 | Additive 2 |
|------------|-------------------|-------------------------|---------------------|------------------------|-----|-------------------------|------------------------|
| A 1 | 2.8% v/v Glycerol | 32% w/v PEG 8000 | None | 100 mM Tris-HCl | 8.5 | 200 mM Cesium Chloride | None |
| A 2 | 5.0% v/v Glycerol | None | None | 100 mM Sodium Citrate | 5.6 | 3.2 M Sodium Formate | None |
| A 3 | 5.0% v/v Glycerol | 20% w/v PEG 4000 | None | None | | None | None |
| A 4 | 10% v/v Glycerol | 15% w/v PEG 5000 MME | 10% w/v Xylitol | None | | None | None |
| A 5 | 10% v/v Glycerol | 28% w/v PEG 8000 | None | 50 mM MES Sodium Salt | 6.5 | 100 mM Ammonium Acetate | None |
| A 6 | 10% v/v Glycerol | 25% w/v PEG 4000 | 4.5% v/v 2-Propanol | 18 mM MES Sodium Salt | 6.5 | None | None |
| B 1 | 10% v/v Glycerol | 20% v/v PEG 400 | 10% w/v PEG 8000 | 50 mM Tris-HCl | 8.5 | 100 mM Sodium Chloride | None |
| B 2 | 10% v/v Glycerol | 15% w/v PEG 6000 | None | 100 mM MES Sodium Salt | 6.5 | 1.0 M Lithium Chloride | None |
| B 3 | 15% v/v Glycerol | 15% w/v PEG 5000 MME | None | None | | None | None |
| B 4 | 15% v/v Glycerol | 20% w/v PEG 6000 | None | 100 mM MES Sodium Salt | 6.5 | None | None |
| B 5 | 15% v/v Glycerol | 30% w/v PEG 8000 | None | 100 mM MES Sodium Salt | 6.5 | 200 mM Sodium Acetate | None |
| B 6 | 15% v/v Glycerol | None | None | 100 mM Tris-HCl | 8.5 | 2.5 M Sodium Chloride | 200 mM Lithium Sulfate |
| C 1 | 15% v/v Glycerol | 15% w/v PEG 20000 | None | None | | None | None |
| C 2 | 15% v/v Glycerol | 35% w/v PEG 4000 | None | 100 mM Tris-Acetate | 8.0 | 100 mM Sodium Citrate | None |
| C 3 | 15% v/v Glycerol | 15% w/v PEG 8000 | None | None | | 500 mM Lithium Sulfate | None |
| C 4 | 20% v/v Glycerol | 16% v/v Ethylene Glycol | 20% w/v PEG 6000 | 30 mM Tris-HCl | 8.5 | 100 mM Sodium Acetate | None |
| C 5 | 20% v/v Glycerol | None | None | None | | 2.0 M Ammonium Sulfate | None |
| C 6 | 20% v/v Glycerol | 20% w/v PEG 4000 | 10% v/v 2-Propanol | 50 mM Sodium Acetate | 4.6 | 100 mM Sodium Chloride | None |
| D 1 | 20% v/v Glycerol | None | None | 100 mM MES Sodium Salt | 6.5 | 2.5 M Ammonium Sulfate | None |
| D 2 | 20% v/v Glycerol | 15% w/v PEG 8000 | None | 100 mM MES Sodium Salt | 6.5 | 160 mM Zinc Acetate | None |
| D 3 | 20% v/v Glycerol | 15% w/v PEG 8000 | None | 80 mM Tris-HCl | 8.5 | 920 mM Sodium Chloride | None |
| D 4 | 20% v/v Glycerol | 6% v/v 2-Propanol | None | None | | 1.8 M Ammonium Sulfate | None |
| D 5 | 20% v/v Glycerol | None | None | 100 mM MES Sodium Salt | 6.5 | 1.8 M Ammonium Sulfate | None |
| D 6 | 20% v/v Glycerol | 20% w/v PEG 6000 | None | 100 mM MES Sodium Salt | 6.5 | None | None |