

## JBScreen Cryo 4

Cat.-No: CC-106

Number	Cryoprotectant 1	Cryoprotectant 2	Cryoprotectant 3	Buffer	pH	Additive 1	Additive 2
<b>A 1</b>	20% v/v Ethylene Glycol	None	None	100 mM HEPES Sodium Salt	7.5	1.6 M Lithium Sulfate	None
<b>A 2</b>	25% v/v Ethylene Glycol	None	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium / Potassium Phosphate	None
<b>A 3</b>	25% v/v Ethylene Glycol	None	None	100 mM HEPES Sodium Salt	7.5	1.0 M Sodium / Potassium Tartrate	None
<b>A 4</b>	40% v/v Ethylene Glycol	None	None	100 mM Sodium Acetate	4.6	None	None
<b>A 5</b>	40% v/v Ethylene Glycol	None	None	100 mM Tris-HCl	8.5	None	None
<b>A 6</b>	40% v/v Ethylene Glycol	None	None	100 mM Imidazole	9.0	200 mM Calcium Acetate	None
<b>B 1</b>	40% v/v Ethylene Glycol	None	None	100 mM MES Sodium Salt	6.5	200 mM Zinc Acetate	None
<b>B 2</b>	50% v/v Ethylene Glycol	None	None	100 mM Tris-HCl	8.5	200 mM Magnesium Chloride	None
<b>B 3</b>	50% v/v Ethylene Glycol	None	None	100 mM HEPES Sodium Salt	7.5	200 mM Lithium Sulfate	None
<b>B 4</b>	35% v/v 2-Propanol	None	None	100 mM Tris-HCl	8.5	None	None
<b>B 5</b>	35% v/v 2-Propanol	None	None	100 mM Sodium Acetate	4.6	None	None
<b>B 6</b>	35% v/v 2-Propanol	None	None	100 mM Imidazole	9.0	200 mM Zinc Acetate	None
<b>C 1</b>	40% v/v 1,2-Propanediol	None	None	100 mM Imidazole	9.0	None	None
<b>C 2</b>	40% v/v 1,2-Propanediol	None	None	100 mM Sodium Acetate	4.6	None	None
<b>C 3</b>	40% v/v 1,2-Propanediol	None	None	100 mM Sodium Citrate	5.6	200 mM Sodium Chloride	None
<b>C 4</b>	40% v/v 1,2-Propanediol	None	None	100 mM Sodium Acetate	4.6	50 mM Calcium Acetate	None
<b>C 5</b>	35% v/v MPD	None	None	100 mM Tris-HCl	8.5	None	None
<b>C 6</b>	30% v/v MPD	None	None	100 mM Sodium Acetate	4.6	20 mM Calcium Chloride	None
<b>D 1</b>	30% v/v MPD	None	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Citrate	None
<b>D 2</b>	30% v/v MPD	None	None	100 mM Sodium Citrate	5.6	200 mM Ammonium Acetate	None
<b>D 3</b>	35% v/v MPD	None	None	100 mM MES Sodium Salt	6.5	200 mM Ammonium Sulfate	None
<b>D 4</b>	40% v/v MPD	None	None	100 mM Sodium Citrate	5.6	None	None
<b>D 5</b>	40% v/v MPD	None	None	100 mM MES Sodium Salt	6.5	200 mM Ammonium Sulfate	None
<b>D 6</b>	40% v/v MPD	None	None	100 mM Imidazole	9.0	200 mM Magnesium Chloride	None