

# Data Sheet

## JBScreen Cryo 3

Cat.-No: CC-105

Number	Cryoprotectant 1	Cryoprotectant 2	Cryoprotectant 3	Buffer	pH	Additive 1	Additive 2
<b>A1</b>	10% v/v PEG 200	24% w/v PEG 3000	None	100 mM Sodium Citrate	5.6	None	None
<b>A2</b>	20% v/v PEG 200	20% w/v PEG 4000	10% v/v 2-Propanol	100 mM HEPES Sodium Salt	7.5	None	None
<b>A3</b>	30% v/v PEG 200	5% w/v PEG 3000	None	100 mM MES Sodium Salt	6.5	None	None
<b>A4</b>	40% v/v PEG 300	5% w/v PEG 1000	None	100 mM Tris-HCl	8.5	None	None
<b>A5</b>	20% v/v PEG 400	30% w/v PEG 4000	None	100 mM Sodium Citrate	5.6	50 mM Ammonium Acetate	None
<b>A6</b>	30% v/v PEG 400	11% w/v PEG 20000	None	100 mM Potassium Phosphate	8.0	None	None
<b>B1</b>	10% v/v PEG 400	20% w/v PEG 5000 MME	None	None		None	None
<b>B2</b>	34% v/v PEG 400	None	None	100 mM HEPES Sodium Salt	7.5	100 mM Sodium Chloride	200 mM Calcium Chloride
<b>B3</b>	40% v/v PEG 400	5% w/v PEG 3000	None	100 mM MES Sodium Salt	6.5	None	None
<b>B4</b>	35% v/v PEG 600	None	None	100 mM PIPES	7.0	100 mM Sodium Chloride	None
<b>B5</b>	30% w/v PEG 1000	None	None	100 mM Phosphate / Citrate	4.4	110 mM Lithium Sulfate	None
<b>B6</b>	40% w/v PEG 1500	None	None	100 mM ADA	6.0	None	None
<b>C 1</b>	17% w/v PEG 1500	10% v/v MPD	None	50 mM PIPES	7.0	None	None
<b>C 2</b>	35% w/v PEG 3350	None	None	None		300 mM Potassium Thiocyanate	None
<b>C 3</b>	25% w/v PEG 4000	18% v/v MPD	None	100 mM Sodium Acetate	4.6	200 mM Ammonium Sulfate	None
<b>C 4</b>	25% w/v PEG 4000	30% v/v Ethylene Glycol	None	None		None	None
<b>C 5</b>	20% w/v PEG 4000	20% v/v MPD	10% v/v 2-Propanol	100 mM Sodium Citrate	5.6	None	None
<b>C 6</b>	10% w/v PEG 4000	20% v/v Ethylene Glycol	None	55 mM ADA	6.0	105 mM Sodium Chloride	None
<b>D 1</b>	2% w/v PEG 4000	25% v/v MPD	None	100 mM HEPES Sodium Salt	7.5	None	None
<b>D 2</b>	20% w/v PEG 8000	20% v/v MPD	None	100 mM HEPES Sodium Salt	7.5	200 mM Rubidium Bromide	None
<b>D 3</b>	28% w/v PEG 8000	None	None	100 mM Glycine	9.5	10 mM Calcium Chloride	None
<b>D 4</b>	10% w/v PEG 8000	15% v/v MPD	None	100 mM MES Sodium Salt	6.5	550 mM Ammonium Sulfate	None
<b>D 5</b>	25% w/v PEG 8000	30% v/v Ethylene Glycol	None	150 mM ADA	6.0	80 mM Ammonium Sulfate	None
<b>D 6</b>	10% w/v PEG 20000	5.6% v/v Ethylene Glycol	30% v/v 2,3-Butanediol	100 mM HEPES Sodium Salt	7.5	None	None