

JBScreen Cryo HTS

Cat.-No.: CC-201S (1.0 ml per well)
CC-201L (1.7 ml per well)

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
A 7	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
A 8	10% v/v Glycerol	15% w/v PEG 6000	None	100 mM MES Sodium Salt	6.5	1.0 M Lithium Chloride	None
A 9	15% v/v Glycerol	15% w/v PEG 5000 MME	None	None		None	None
A 10	15% v/v Glycerol	20% w/v PEG 6000	None	100 mM MES Sodium Salt	6.5	None	None
A 11	15% v/v Glycerol	30% w/v PEG 8000	None	100 mM MES Sodium Salt	6.5	200 mM Sodium Acetate	None
A 12	15% v/v Glycerol	None	None	100 mM Tris-HCl	8.5	2.5 M Sodium Chloride	200 mM Lithium Sulfate
B1	15% v/v Glycerol	15% w/v PEG 20000	None	None		None	None
B2	15% v/v Glycerol	35% w/v PEG 4000	None	100 mM Tris-Acetate	8.0	200 mM Sodium Citrate	None
B3	15% v/v Glycerol	15% w/v PEG 8000	None	None		500 mM Lithium Sulfate	None
B4	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
B5	20% v/v Glycerol	None	None	None		2.0 M Ammonium Sulfate	None
B6	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
B7	20% v/v Glycerol	None	None	100 mM MES Sodium Salt	6.5	2.5 M Ammonium Sulfate	None
B8	20% v/v Glycerol	15% w/v PEG 8000	None	100 mM MES Sodium Salt	6.5	160 mM Zinc Acetate	None
B9	20% v/v Glycerol	15% w/v PEG 8000	None	80 mM Tris-HCl	8.5	920 mM Sodium Chloride	None
B10	20% v/v Glycerol	6% v/v 2-Propanol	None	None		1.8 M Ammonium Sulfate	None
B11	20% v/v Glycerol	None	None	100 mM MES Sodium Salt	6.5	1.8 M Ammonium Sulfate	None
B12	20% v/v Glycerol	20% w/v PEG 6000	None	100 mM MES Sodium Salt	6.5	None	None

Data Sheet

JBScreen Cryo HTS

Cat.-No.: CC-201S (1.0 ml per well)

CC-201L (1.7 ml per well)

Number	Cryoprotectant 1	Cryoprotectant 2	Cryoprotectant 3	Buffer	pH	Additive 1	Additive 2
C 1	20% v/v Glycerol	16% w/v PEG 8000	None	80 mM MES Sodium Salt	6.5	160 mM Magnesium Acetate	None
C 2	20% v/v Glycerol	7% w/v PEG 8000	None	100 mM MES Sodium Salt	6.5	None	None
C 3	24% v/v Glycerol	None	None	None		2.3 M Ammonium Sulfate	None
C 4	25% v/v Glycerol	None	None	None		750 mM Sodium Bromide	None
C 5	25% v/v Glycerol	None	None	100 mM Sodium Citrate	5.6	1.2 M Ammonium dihydrogen Phosphate	None
C 6	25% v/v Glycerol	None	None	None		1.5 M Ammonium Sulfate	None
C 7	25% v/v Glycerol	None	None	100 mM MES Sodium Salt	6.5	2.2 M Ammonium Sulfate	None
C 8	25% v/v Glycerol	None	None	100 mM Sodium / Potassium Phosphate	6.5	2.5 M Ammonium Sulfate	None
C 9	30% v/v Glycerol	None	None	None	7.0	1.26 M Ammonium dihydrogen Phosphate	None
C 10	30% v/v Glycerol	None	None	100 mM Sodium Acetate	4.6	2.0 M Ammonium Sulfate	None
C 11	30% v/v Glycerol	None	None	50 mM MES Sodium Salt	6.5	2.3 M Ammonium Sulfate	None
C 12	30% v/v Glycerol	None	None	None		2.0 M Ammonium Sulfate	None
D 1	30% v/v Glycerol	15% w/v PEG 4000	5% v/v 2-Propanol	100 mM HEPES Sodium Salt	7.5	None	None
D 2	30% v/v Glycerol	20% w/v PEG 4000	20% v/v 2-Propanol	100 mM Tris-HCl	8.5	None	None
D 3	30% v/v Glycerol	None	None	50 mM MES Sodium Salt	6.5	1.2 M Ammonium Sulfate	None
D 4	30% v/v Glycerol	15% w/v PEG 6000	None	500 mM Tris-HCl	8.5	None	None
D 5	30% v/v Glycerol	None	None	None		2.0 M Ammonium Sulfate	None
D 6	30% v/v Glycerol	30% w/v PEG 1500	None	50 mM MES Sodium Salt	6.5	None	None
D 7	30% v/v Glycerol	None	None	100 mM Sodium / Potassium Phosphate	6.5	1.0 M Ammonium Sulfate	None
D 8	30% v/v Glycerol	30% w/v PEG 4000	None	100 mM Sodium Acetate	4.6	None	None
D 9	30% v/v Glycerol	None	None	100 mM MES Sodium Salt	6.5	1.4 M Sodium Acetate	None
D 10	30% v/v Glycerol	None	None	150 mM HEPES Sodium Salt	7.5	1.6 M Sodium dihydrogen Phosphate	450 mM di-Potassium hydrogen Phosphate
D 11	35% v/v Glycerol	None	None	20 mM HEPES Sodium Salt	7.5	None	None
D 12	35% v/v Glycerol	7% w/v PEG 6000	None	50 mM MES Sodium Salt	6.5	None	None

JBScreen Cryo HTS

Cat.-No.: CC-201S (1.0 ml per well)
CC-201L (1.7 ml per well)

Number	Cryoprotectant 1	Cryoprotectant 2	Cryoprotectant 3	Buffer	pH	Additive 1	Additive 2
E1	10% v/v PEG 200	24% w/v PEG 3000	None	100 mM Sodium Citrate	5.6	None	None
E2	20% v/v PEG 200	20% w/v PEG 4000	10% v/v 2-Propanol	100 mM HEPES Sodium Salt	7.5	None	None
E3	30% v/v PEG 200	5% w/v PEG 3000	None	100 mM MES Sodium Salt	6.5	None	None
E4	40% v/v PEG 300	5% w/v PEG 1000	None	100 mM Tris-HCl	8.5	None	None
E5	20% v/v PEG 400	30% w/v PEG 4000	None	100 mM Sodium Citrate	5.6	50 mM Ammonium Acetate	None
E6	30% v/v PEG 400	11% w/v PEG 20000	None	100 mM Potassium hydrogen Phosphate	8.0	None	None
E7	10% v/v PEG 400	20% w/v PEG 5000 MME	None	None		None	None
E8	34% v/v PEG 400	None	None	100 mM HEPES Sodium Salt	7.5	100 mM Sodium Chloride	200 mM Calcium Chloride
E9	40% v/v PEG 400	5% w/v PEG-3000	None	100 mM MES Sodium Salt	6.5	None	None
E10	35% v/v PEG 600	None	None	100 mM PIPES	7.0	100 mM Sodium Chloride	None
E11	30% w/v PEG 1000	None	None	100 mM Phosphate / Citrate	4.4	110 mM Lithium Sulfate	None
E12	40% w/v PEG 1500	None	None	100 mM ADA	6.0	None	None
F 1	17% w/v PEG 1500	10% v/v MPD	None	50 mM PIPES	7.0	None	None
F 2	35% w/v PEG 3350	None	None	None		300 mM Potassium Thiocyanate	None
F 3	25% w/v PEG 4000	18% v/v MPD	None	100 mM Sodium Acetate	4.6	200 mM Ammonium Sulfate	None
F 4	25% w/v PEG 4000	30% v/v Ethylene Glycol	None	None		None	None
F 5	20% w/v PEG 4000	20% v/v MPD	10% v/v 2-Propanol	100 mM Sodium Citrate	5.6	None	None
F 6	10% w/v PEG 4000	20% v/v Ethylene Glycol	None	55 mM ADA	6.0	105 mM Sodium Chloride	None
F 7	2% w/v PEG 4000	25% v/v MPD	None	100 mM HEPES Sodium Salt	7.5	None	None
F 8	20% w/v PEG 8000	20% v/v MPD	None	100 mM HEPES Sodium Salt	7.5	200 mM Rubidium Bromide	None
F 9	28% w/v PEG 8000	None	None	100 mM Glycine	9.5	10 mM Calcium Chloride	None
F 10	10% w/v PEG 8000	15% v/v MPD	None	100 mM MES Sodium Salt	6.5	550 mM Ammonium Sulfate	None
F 11	25 %/v PEG 8000	30% v/v Ethylene Glycol	None	150 mM ADA	6.0	80 mM Ammonium Sulfate	None
F 12	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

JBScreen Cryo HTS

Cat.-No.: CC-201S (1.0 ml per well)
CC-201L (1.7 ml per well)

Number	Cryoprotectant 1	Cryoprotectant 2	Cryoprotectant 3	Buffer	pH	Additive 1	Additive 2
G 1	20% v/v Ethylene Glycol	None	None	100 mM HEPES Sodium Salt	7.5	1.6 M Lithium Sulfate	None
G 2	25% v/v Ethylene Glycol	None	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium / Potassium Phosphate	None
G 3	25% v/v Ethylene Glycol	None	None	100 mM HEPES Sodium Salt	7.5	1.0 M Sodium / Potassium	None
G 4	40% v/v Ethylene Glycol	None	None	100 mM Sodium Acetate	4.6	None	None
G 5	40% v/v Ethylene Glycol	None	None	100 mM Tris-HCl	8.5	None	None
G 6	40% v/v Ethylene Glycol	None	None	100 mM Imidazole	9.0	200 mM Calcium Acetate	None
G 7	40% v/v Ethylene Glycol	None	None	100 mM MES Sodium Salt	6.5	200 mM Zinc Acetate	None
G 8	50% v/v Ethylene Glycol	None	None	100 mM Tris-HCl	8.5	200 mM Magnesium Chloride	None
G 9	50% v/v Ethylene Glycol	None	None	100 mM HEPES Sodium Salt	7.5	200 mM Lithium Sulfate	None
G 10	35% v/v 2-Propanol	None	None	100 mM Tris-HCl	8.5	None	None
G 11	35% v/v 2-Propanol	None	None	100 mM Sodium Acetate	4.6	None	None
G 12	35% v/v 2-Propanol	None	None	100 mM Imidazole	9.0	200 mM Zinc Acetate	None
H 1	40% v/v 1,2-Propanediol	None	None	100 mM Imidazole	9.0	None	None
H 2	40% v/v 1,2-Propanediol	None	None	100 mM Sodium Acetate	4.6	None	None
H 3	40% v/v 1,2-Propanediol	None	None	100 mM Sodium Citrate	5.6	200 mM Sodium Chloride	None
H 4	40% v/v 1,2-Propanediol	None	None	100 mM Sodium Acetate	4.6	50 mM Calcium Acetate	None
H 5	35% v/v MPD	None	None	100 mM Tris-HCl	8.5	None	None
H 6	30% v/v MPD	None	None	100 mM Sodium Acetate	4.6	20 mM Calcium Chloride	None
H 7	30% v/v MPD	None	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Citrate	None
H 8	30% v/v MPD	None	None	100 mM Sodium Citrate	5.6	200 mM Ammonium Acetate	None
H 9	35% v/v MPD	None	None	100 mM MES Sodium Salt	6.5	200 mM Ammonium Sulfate	None
H 10	40% v/v MPD	None	None	100 mM Sodium Citrate	5.6	None	None
H 11	40% v/v MPD	None	None	100 mM MES Sodium Salt	6.5	200 mM Ammonium Sulfate	None
H 12	40% v/v MPD	None	None	100 mM Imidazole	9.0	200 mM Magnesium Chloride	None