

## JBScreen Classic HTS II

Cat.-No.: CS-202S (1.0 ml per well)

CS-202L (1.7 ml per well)

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
<b>A 1</b>	500 mM Ammonium Sulfate	1.0 M Lithium Sulfate	None		100 mM Sodium Citrate
<b>A 2</b>	1.0 M Ammonium Sulfate	None	100 mM Sodium Acetate	4.6	None
<b>A 3</b>	1.0 M Ammonium Sulfate	2 % w/v PEG 400	100 mM HEPES Sodium Salt	7.5	None
<b>A 4</b>	1.0 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	None
<b>A 5</b>	1.2 M Ammonium Sulfate	3 % w/v 2-Propanol	None		50 mM Sodium Citrate
<b>A 6</b>	1.5 M Ammonium Sulfate	15 % w/v Glycerol	100 mM Tris-HCl	8.5	None
<b>A 7</b>	1.6 M Ammonium Sulfate	1.0 M Lithium Sulfate	None		None
<b>A 8</b>	1.6 M Ammonium Sulfate	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Chloride
<b>A 9</b>	1.6 M Ammonium Sulfate	2 % w/v PEG 1000	100 mM HEPES Sodium Salt	7.5	None
<b>A 10</b>	1.8 M Ammonium Sulfate	None	100 mM MES Sodium Salt	6.5	None
<b>A 11</b>	2.0 M Ammonium Sulfate	2.0 M Sodium Chloride	None		None
<b>A 12</b>	2.0 M Ammonium Sulfate	None	100 mM Sodium Acetate	4.6	None
<b>B 1</b>	2.0 M Ammonium Sulfate	5 % w/v PEG 400	100 mM MES Sodium Salt	6.5	None
<b>B 2</b>	2.0 M Ammonium Sulfate	None	100 mM Tris-HCl	8.5	None
<b>B 3</b>	2.2 M Ammonium Sulfate	20 % w/v Glycerol	None		None
<b>B 4</b>	2.4 M Ammonium Sulfate	None			100 mM Sodium Citrate
<b>B 5</b>	3.0 M Ammonium Sulfate	1 % w/v MPD	None		None
<b>B 6</b>	3.0 M Ammonium Sulfate	10 % w/v Glycerol	None		None
<b>B 7</b>	3.5 M Ammonium Sulfate	None	100 mM HEPES Sodium Salt	7.5	None
<b>B 8</b>	3.5 M Ammonium Sulfate	1 % w/v MPD	100 mM MES Sodium Salt	6.5	None
<b>B 9</b>	10 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	100 mM Sodium Citrate
<b>B 10</b>	12 % w/v MPD	None	100 mM Tris-HCl	8.5	50 mM Magnesium Chloride
<b>B 11</b>	15 % w/v MPD	5 % w/v PEG 4000	100 mM Imidazole-HCl	8.0	None
<b>B 12</b>	15 % w/v MPD	None	100 mM Sodium Citrate	5.6	200 mM Ammonium Acetate

# Data Sheet

## JBScreen Classic HTS II

Cat.-No.: CS-202S (1.0 ml per well)

CS-202L (1.7 ml per well)

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
<b>C 1</b>	15 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Citrate	None
<b>C 2</b>	20 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	100 mM Sodium Citrate	None
<b>C 3</b>	20 % w/v MPD	None	100 mM Imidazole-HCl	8.0	None	None
<b>C 4</b>	20 % w/v MPD	4 % w/v Glycerol	None		200 mM Sodium Chloride	None
<b>C 5</b>	30 % w/v MPD	None	100 mM Sodium Acetate	4.6	20 mM Calcium Chloride	None
<b>C 6</b>	30 % w/v MPD	None	100 mM Sodium Citrate	5.6	200 mM Ammonium Acetate	None
<b>C 7</b>	30 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	500 mM Ammonium Sulfate	None
<b>C 8</b>	30 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Citrate	None
<b>C 9</b>	30 % w/v MPD	10 % w/v PEG 4000	100 mM Imidazole-HCl	8.0	None	None
<b>C 10</b>	30 % w/v MPD	20 % w/v Ethanol	None		None	None
<b>C 11</b>	35 % w/v MPD	None	100 mM Imidazole-HCl	8.0	None	None
<b>C 12</b>	40 % w/v MPD	None	100 mM Tris-HCl	8.5	None	None
<b>D 1</b>	47 % w/v MPD	None	100 mM HEPES Sodium Salt	7.5	None	None
<b>D 2</b>	47 % w/v MPD	2 % w/v tert.-Butanol	None		None	None
<b>D 3</b>	50 % w/v MPD	None	None		None	None
<b>D 4</b>	50 % w/v MPD	20 % w/v 2-Propanol	None		50 mM Sodium Chloride	50 mM Sodium Acetate
<b>D 5</b>	50 % w/v MPD	None	100 mM Tris-HCl	8.5	100 mM Ammonium dihydrogen Phosphate	None
<b>D 6</b>	55 % w/v MPD	None	None		None	None
<b>D 7</b>	60 % w/v MPD	None	100 mM Sodium Acetate	4.6	10 mM Calcium Chloride	None
<b>D 8</b>	70 % w/v MPD	None	100 mM MES Sodium Salt	6.5	None	None
<b>D 9</b>	70 % w/v MPD	None	100 mM Tris-HCl	8.5	None	None
<b>D 10</b>	20 % w/v Methanol	None	100 mM Tris-HCl	8.5	10 mM Calcium Chloride	None
<b>D 11</b>	2 % w/v Ethanol	None	100 mM Tris-HCl	8.5	None	None
<b>D 12</b>	5 % w/v Ethanol	5 % w/v MPD	100 mM Tris-HCl	8.5	200 mM Sodium Chloride	None

## JBScreen Classic HTS II

Cat.-No.: CS-202S (1.0 ml per well)

CS-202L (1.7 ml per well)

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
<b>E 1</b>	10 % w/v Ethanol	None	100 mM Tris-HCl	8.5	None
<b>E 2</b>	12 % w/v Ethanol	4 % w/v PEG 400	100 mM Sodium Acetate	4.6	None
<b>E 3</b>	14 % w/v Ethanol	5 % w/v Glycerol	100 mM Tris-HCl	8.5	None
<b>E 4</b>	18 % w/v Ethanol	None	100 mM Tris-HCl	8.5	None
<b>E 5</b>	20 % w/v Ethanol	10 % w/v Glycerol	None		None
<b>E 6</b>	30 % w/v Ethanol	10 % w/v PEG 6000	None		100 mM Sodium Acetate
<b>E 7</b>	45 % w/v Ethanol	None	None		None
<b>E 8</b>	50 % w/v Ethanol	None	None		10 mM Sodium Acetate
<b>E 9</b>	60 % w/v Ethanol	1.5 % w/v PEG 6000	None		50 mM Sodium Acetate
<b>E 10</b>	60 % w/v Ethanol	None	None		100 mM Sodium Chloride
<b>E 11</b>	5 %wV 2-Propanol	None	100 mM HEPES Sodium Salt	7.5	
<b>E 12</b>	10 % w/v 2-Propanol	None	100 mM Sodium Acetate	4.6	200 mM Calcium Chloride
<b>F 1</b>	10 % w/v 2-Propanol	None	100 mM HEPES Sodium Salt	7.5	200 mM Sodium Citrate
<b>F 2</b>	10 % w/v 2-Propanol	None	100 mM Tris-HCl	8.5	10 mM Magnesium Chloride
<b>F 3</b>	15 % w/v 2-Propanol	None	100 mM MES Sodium Salt	6.5	200 mM Sodium Citrate
<b>F 4</b>	15 % w/v 2-Propanol	None	100 mM HEPES Sodium Salt	7.5	200 mM Magnesium Chloride
<b>F 5</b>	15 % w/v 2-Propanol	None	100 mM Tris-HCl	8.5	200 mM Ammonium Acetate
<b>F 6</b>	20 % w/v 2-Propanol	None	100 mM Sodium Acetate	4.6	200 mM Calcium Chloride
<b>F 7</b>	25 % w/v 2-Propanol	None	100 mM HEPES Sodium Salt	7.5	100 mM Magnesium Chloride
<b>F 8</b>	30 % w/v 2-Propanol	None	100 mM MES Sodium Salt	6.5	200 mM Sodium Citrate
<b>F 9</b>	30 % w/v 2-Propanol	None	100 mM Tris-HCl	8.5	200 mM Ammonium Acetate
<b>F 10</b>	25 % w/v tert-Butanol	None	100 mM Tris-HCl	8.5	100 mM Calcium Chloride
<b>F 11</b>	35 % w/v tert-Butanol	None	100 mM Sodium Citrate	5.6	None
<b>F 12</b>	200 mM Ammonium dihydrogen Phosphate	None	None		None

## JBScreen Classic HTS II

Cat.-No.: CS-202S (1.0 ml per well)

CS-202L (1.7 ml per well)

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive
<b>G 1</b>	200 mM Potassium / Sodium Tartrate	None	None		None
<b>G 2</b>	200 mM Magnesium Acetate	None	None		None
<b>G 3</b>	400 mM Ammonium dihydrogen Phosphate	None	None		None
<b>G 4</b>	400 mM Potassium / Sodium Tartrate	None	100 mM Tris-HCl	8.5	None
<b>G 5</b>	500 mM Ammonium dihydrogen Phosphate	None	None		200 mM Sodium Citrate
<b>G 6</b>	500 mM Sodium Acetate	None	100 mM Imidazole-HCl	8.0	None
<b>G 7</b>	700 mM Sodium Citrate	None	100 mM HEPES Sodium Salt	7.5	None
<b>G 8</b>	800 mM Potassium / Sodium Tartrate	None	100 mM HEPES Sodium Salt	7.5	None
<b>G 9</b>	1.0 M Ammonium dihydrogen Phosphate	None	100 mM Sodium Citrate	5.6	None
<b>G 10</b>	1.0 M Ammonium dihydrogen Phosphate	None	100 mM Tris-HCl	8.5	None
<b>G 11</b>	1.0 M Lithium Sulfate	None	100 mM Tris-HCl	8.5	10 mM Nickel (II) Chloride
<b>G 12</b>	1.0 M Sodium Acetate	None	100 mM Imidazole-HCl	8.0	None
<b>H 1</b>	1.4 M Sodium Acetate	None	100 mM MES Sodium Salt	6.5	None
<b>H 2</b>	1.5 M Lithium Sulfate	None	100 mM Tris-HCl	8.5	None
<b>H 3</b>	1.5 M Sodium Citrate	None	None	6.5	None
<b>H 4</b>	1.6 M Magnesium Sulfate	None	100 mM MES Sodium Salt	6.5	None
<b>H 5</b>	1.6 M Potassium / Sodium Tartrate	None	100 mM MES Sodium Salt	6.5	None
<b>H 6</b>	2.0 M Ammonium Formate	None	100 mM MES Sodium Salt	6.5	None
<b>H 7</b>	2.0 M Ammonium dihydrogen Phosphate	None	100 mM Tris-HCl	8.5	None
<b>H 8</b>	2.0 M Sodium Chloride	None	100 mM MES Sodium Salt	6.5	200 mM Sodium Acetate
<b>H 9</b>	2.0 M Sodium Formate	None	100 mM Sodium Acetate	4.6	None
<b>H 10</b>	1.0 M Ammonium dihydrogen Phosphate	30 % w/v Glycerol	100 mM Tris-HCl	8.5	None
<b>H 11</b>	4.0 M Sodium Chloride	None	100 mM HEPES Sodium Salt	7.5	None
<b>H 12</b>	3.0 M Sodium Formate	None	None		None