

## JBScreen PEG/Salt HTS

Cat.-No.: CS-205S (1.0 ml per well)

CS-205L (1.7 ml per well)

Number	Precipitant	Salt
<b>A 1</b>	20% w/v PEG 3350	200 mM Ammonium Acetate
<b>A 2</b>	20% w/v PEG 3350	200 mM Ammonium Chloride
<b>A 3</b>	20% w/v PEG 3350	200 mM Ammonium Fluoride
<b>A 4</b>	20% w/v PEG 3350	200 mM Ammonium Formate
<b>A 5</b>	20% w/v PEG 3350	200 mM Ammonium Iodide
<b>A 6</b>	20% w/v PEG 3350	200 mM Ammonium Nitrate
<b>A 7</b>	20% w/v PEG 3350	200 mM Ammonium dihydrogen Phosphate
<b>A 8</b>	20% w/v PEG 3350	200 mM di-Ammonium hydrogen Phosphate
<b>A 9</b>	20% w/v PEG 3350	200 mM Ammonium Sulfate
<b>A 10</b>	20% w/v PEG 3350	200 mM Ammonium Sulfite
<b>A 11</b>	20% w/v PEG 3350	200 mM Calcium Acetate
<b>A 12</b>	20% w/v PEG 3350	200 mM Calcium Chloride
<b>B 1</b>	20% w/v PEG 3350	200 mM di-Ammonium Tartrate
<b>B 2</b>	20% w/v PEG 3350	200 mM Potassium Formate
<b>B 3</b>	20% w/v PEG 3350	200 mM Lithium Acetate
<b>B 4</b>	20% w/v PEG 3350	200 mM Lithium Chloride
<b>B 5</b>	20% w/v PEG 3350	200 mM Lithium Citrate
<b>B 6</b>	20% w/v PEG 3350	200 mM Lithium Nitrate
<b>B 7</b>	20% w/v PEG 3350	200 mM Lithium Sulfate
<b>B 8</b>	20% w/v PEG 3350	200 mM Magnesium Acetate
<b>B 9</b>	20% w/v PEG 3350	200 mM Magnesium Chloride
<b>B 10</b>	20% w/v PEG 3350	200 mM Magnesium Formate
<b>B 11</b>	20% w/v PEG 3350	200 mM Magnesium Nitrate
<b>B 12</b>	20% w/v PEG 3350	200 mM Magnesium Sulfate

## JBScreen PEG/Salt HTS

Cat.-No.: CS-205S (1.0 ml per well)

CS-205L (1.7 ml per well)

Number	Precipitant	Salt
<b>C 1</b>	20% w/v PEG 3350	200 mM Potassium Acetate
<b>C 2</b>	20% w/v PEG 3350	200 mM Potassium Chloride
<b>C 3</b>	20% w/v PEG 3350	200 mM Potassium Fluoride
<b>C 4</b>	20% w/v PEG 3350	200 mM Potassium Iodide
<b>C 5</b>	20% w/v PEG 3350	200 mM Potassium Nitrate
<b>C 6</b>	20% w/v PEG 3350	200 mM Potassium dihydrogen Phosphate
<b>C 7</b>	20% w/v PEG 3350	200 mM di-Potassium hydrogen Phosphate
<b>C 8</b>	20% w/v PEG 3350	200 mM Potassium Sulfate
<b>C 9</b>	20% w/v PEG 3350	200 mM Potassium Thiocyanate
<b>C 10</b>	20% w/v PEG 3350	200 mM Potassium / Sodium Tartrate
<b>C 11</b>	20% w/v PEG 3350	200 mM Sodium Acetate
<b>C 12</b>	20% w/v PEG 3350	200 mM Sodium Chloride
<b>D 1</b>	20% w/v PEG 3350	200 mM Sodium Citrate
<b>D 2</b>	20% w/v PEG 3350	200 mM Sodium Fluoride
<b>D 3</b>	20% w/v PEG 3350	200 mM Sodium Formate
<b>D 4</b>	20% w/v PEG 3350	200 mM Sodium Iodide
<b>D 5</b>	20% w/v PEG 3350	200 mM Sodium Isothiocyanate
<b>D 6</b>	20% w/v PEG 3350	200 mM Sodium Nitrate
<b>D 7</b>	20% w/v PEG 3350	200 mM Sodium dihydrogen Phosphate
<b>D 8</b>	20% w/v PEG 3350	200 mM di-Sodium hydrogen Phosphate
<b>D 9</b>	20% w/v PEG 3350	200 mM Sodium Sulfate
<b>D 10</b>	20% w/v PEG 3350	200 mM di-Sodium Tartrate
<b>D 11</b>	20% w/v PEG 3350	200 mM tri-Potassium Citrate
<b>D 12</b>	20% w/v PEG 3350	200 mM Zinc Acetate

## JBScreen PEG/Salt HTS

Cat.-No.: CS-205S (1.0 ml per well)

CS-205L (1.7 ml per well)

Number	Precipitant	Salt
<b>E 1</b>	20 % w/v PEG 5000 MME	200 mM Ammonium Acetate
<b>E 2</b>	20 % w/v PEG 5000 MME	200 mM Ammonium Chloride
<b>E 3</b>	20 % w/v PEG 5000 MME	200 mM Ammonium Fluoride
<b>E 4</b>	20 % w/v PEG 5000 MME	200 mM Ammonium Formate
<b>E 5</b>	20 % w/v PEG 5000 MME	200 mM Ammonium Iodide
<b>E 6</b>	20 % w/v PEG 5000 MME	200 mM Ammonium Nitrate
<b>E 7</b>	20 % w/v PEG 5000 MME	200 mM Ammonium dihydrogen Phosphate
<b>E 8</b>	20 % w/v PEG 5000 MME	200 mM di-Ammonium hydrogen Phosphate
<b>E 9</b>	20 % w/v PEG 5000 MME	200 mM Ammonium Sulfate
<b>E 10</b>	20 % w/v PEG 5000 MME	200 mM Ammonium Sulfite
<b>E 11</b>	20 % w/v PEG 5000 MME	200 mM Calcium Acetate
<b>E 12</b>	20 % w/v PEG 5000 MME	200 mM Calcium Chloride
<b>F 1</b>	20 % w/v PEG 5000 MME	200 mM di-Ammonium Tartrate
<b>F 2</b>	20 % w/v PEG 5000 MME	200 mM Potassium Formate
<b>F 3</b>	20 % w/v PEG 5000 MME	200 mM Lithium Acetate
<b>F 4</b>	20 % w/v PEG 5000 MME	200 mM Lithium Chloride
<b>F 5</b>	20 % w/v PEG 5000 MME	200 mM Lithium Citrate
<b>F 6</b>	20 % w/v PEG 5000 MME	200 mM Lithium Nitrate
<b>F 7</b>	20 % w/v PEG 5000 MME	200 mM Lithium Sulfate
<b>F 8</b>	20 % w/v PEG 5000 MME	200 mM Magnesium Acetate
<b>F 9</b>	20 % w/v PEG 5000 MME	200 mM Magnesium Chloride
<b>F 10</b>	20 % w/v PEG 5000 MME	200 mM Magnesium Formate
<b>F 11</b>	20 % w/v PEG 5000 MME	200 mM Magnesium Nitrate
<b>F 12</b>	20 % w/v PEG 5000 MME	200 mM Magnesium Sulfate

## JBScreen PEG/Salt HTS

Cat.-No.: CS-205S (1.0 ml per well)

CS-205L (1.7 ml per well)

Number	Precipitant	Salt
<b>G 1</b>	20 % w/v PEG 5000 MME	200 mM Potassium Acetate
<b>G 2</b>	20 % w/v PEG 5000 MME	200 mM Potassium Chloride
<b>G 3</b>	20 % w/v PEG 5000 MME	200 mM Potassium Fluoride
<b>G 4</b>	20 % w/v PEG 5000 MME	200 mM Potassium Iodide
<b>G 5</b>	20 % w/v PEG 5000 MME	200 mM Potassium Nitrate
<b>G 6</b>	20 % w/v PEG 5000 MME	200 mM Potassium dihydrogen Phosphate
<b>G 7</b>	20 % w/v PEG 5000 MME	200 mM di-Potassium hydrogen Phosphate
<b>G 8</b>	20 % w/v PEG 5000 MME	200 mM Potassium Sulfate
<b>G 9</b>	20 % w/v PEG 5000 MME	200 mM Potassium Thiocyanate
<b>G 10</b>	20 % w/v PEG 5000 MME	200 mM Potassium / Sodium Tartrate
<b>G 11</b>	20 % w/v PEG 5000 MME	200 mM Sodium Acetate
<b>G 12</b>	20 % w/v PEG 5000 MME	200 mM Sodium Chloride
<b>H 1</b>	20 % w/v PEG 5000 MME	200 mM Sodium Citrate
<b>H 2</b>	20 % w/v PEG 5000 MME	200 mM Sodium Fluoride
<b>H 3</b>	20 % w/v PEG 5000 MME	200 mM Sodium Formate
<b>H 4</b>	20 % w/v PEG 5000 MME	200 mM Sodium Iodide
<b>H 5</b>	20 % w/v PEG 5000 MME	200 mM Sodium Isothiocyanate
<b>H 6</b>	20 % w/v PEG 5000 MME	200 mM Sodium Nitrate
<b>H 7</b>	20 % w/v PEG 5000 MME	200 mM Sodium dihydrogen Phosphate
<b>H 8</b>	20 % w/v PEG 5000 MME	200 mM di-Sodium hydrogen Phosphate
<b>H 9</b>	20 % w/v PEG 5000 MME	200 mM Sodium Sulfate
<b>H 10</b>	20 % w/v PEG 5000 MME	200 mM di-Sodium Tartrate
<b>H 11</b>	20 % w/v PEG 5000 MME	200 mM tri-Potassium Citrate
<b>H 12</b>	20 % w/v PEG 5000 MME	200 mM Zinc Acetate