

## JBScreen Classic 8

Cat.-No: CS-108L

Number	Precipitant 1	Precipitant 2	Buffer	pH	Additive 1	Additive 2
<b>A 1</b>	50 % w/v MPD	15 % w/v Ethanol	None		10 mM Sodium Acetate	None
<b>A 2</b>	50 % w/v MPD	20 % w/v 2-Propanol	None		50 mM Sodium Chloride	50 mM Sodium Acetate
<b>A 3</b>	50 % w/v MPD	None	100 mM Tris-HCl	8.5	100 mM Ammonium dihydrogen Phosphate	None
<b>A 4</b>	55 % w/v MPD	None	None		None	None
<b>A 5</b>	60 % w/v MPD	None	100 mM Sodium Acetate	4.6	10 mM Calcium Chloride	None
<b>A 6</b>	60 % w/v MPD	None	None		20 mM Sodium Acetate	None
<b>B 1</b>	70 % w/v MPD	None	100 mM MES Sodium Salt	6.5	None	None
<b>B 2</b>	70 % w/v MPD	None	100 mM Tris-HCl	8.5	None	None
<b>B 3</b>	20 % w/v Methanol	None	100 mM Tris-HCl	8.5	10 mM Calcium Chloride	None
<b>B 4</b>	2 % w/v Ethanol	None	100 mM Tris-HCl	8.5	None	None
<b>B 5</b>	5 % w/v Ethanol	5 % w/v MPD	100 mM HEPES Sodium Salt	7.5	None	None
<b>B 6</b>	5 % w/v Ethanol	5 % w/v MPD	100 mM Tris-HCl	8.5	200 mM Sodium Chloride	None
<b>C 1</b>	10 % w/v Ethanol	None	100 mM Tris-HCl	8.5	None	None
<b>C 2</b>	12 % w/v Ethanol	4 % w/v PEG 400	100 mM Sodium Acetate	4.6	None	None
<b>C 3</b>	14 % w/v Ethanol	5 % w/v Glycerol	100 mM Tris-HCl	8.5	None	None
<b>C 4</b>	18 % w/v Ethanol	None	100 mM Tris-HCl	8.5	None	None
<b>C 5</b>	20 % w/v Ethanol	None	None		None	None
<b>C 6</b>	20 % w/v Ethanol	10 % w/v Glycerol	None		None	None
<b>D 1</b>	30 % w/v Ethanol	10 % w/v PEG 6000	None		100 mM Sodium Acetate	None
<b>D 2</b>	45 % w/v Ethanol	None	None		None	None
<b>D 3</b>	50 % w/v Ethanol	None	None		10 mM Sodium Acetate	None
<b>D 4</b>	60 % w/v Ethanol	1.5 % w/v PEG 6000	None		50 mM Sodium Acetate	None
<b>D 5</b>	60 % w/v Ethanol	None	None		100 mM Sodium Chloride	None
<b>D 6</b>	2 % w/v 2-Propanol	None	100 mM Tris-HCl	8.5	10 mM Magnesium Sulfate	None