

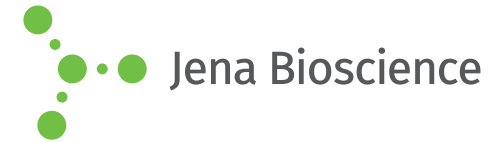


## JBScreen Classic 10

(Salt based)

Cat.-No.: CS-110L

## SCREEN FORMULATION



Jena Bioscience

No.	Precipitant 1	Precipitant 2	Buffer	Additive
10/A1	500 mM Sodium acetate	none	100 mM Imidazole; pH 8.0	none
10/A2	700 mM tri-Sodium citrate	none	100 mM HEPES; pH 7.5	none
10/A3	700 mM Lithium sulfate	none	100 mM TRIS; pH 8.5	none
10/A4	800 mM Potassium Sodium tartrate	none	100 mM HEPES; pH 7.5	none
10/A5	1 M Ammonium di-hydrogen phosphate	none	100 mM tri-Sodium citrate; pH 5.6	none
10/A6	1 M Ammonium di-hydrogen phosphate	none	100 mM TRIS; pH 8.5	none
10/B1	1 M Lithium sulfate	none	100 mM TRIS; pH 8.5	10 mM Nickel (II) chloride
10/B2	1 M Sodium acetate	none	100 mM Imidazole; pH 8.0	none
10/B3	1 M Sodium formate	none	100 mM Sodium acetate; pH 4.6	none
10/B4	1.4 M Sodium acetate	none	100 mM MES; pH 6.5	none
10/B5	1.4 M tri-Sodium citrate	none	100 mM HEPES; pH 7.5	none
10/B6	1.5 M Lithium sulfate	none	100 mM TRIS; pH 8.5	none
10/C1	1.5 M tri-Sodium citrate; pH 6.5	none	none	none
10/C2	1.6 M Magnesium sulfate	none	100 mM MES; pH 6.5	none
10/C3	1.6 M Potassium Sodium tartrate	none	100 mM MES; pH 6.5	none
10/C4	2 M Ammonium formate	none	100 mM MES; pH 6.5	none
10/C5	2 M Ammonium di-hydrogen phosphate	none	100 mM TRIS; pH 8.5	none
10/C6	2 M Sodium formate	none	none	none
10/D1	2 M Magnesium chloride	none	100 mM TRIS; pH 8.5	none
10/D2	2 M Sodium chloride	none	100 mM MES; pH 6.5	200 mM Sodium acetate
10/D3	2 M Sodium formate	none	100 mM Sodium acetate; pH 4.6	none
10/D4	1 M Ammonium di-hydrogen phosphate	30 % w/v Glycerol	100 mM TRIS; pH 8.5	none
10/D5	4 M Sodium chloride	none	100 mM HEPES; pH 7.5	none
10/D6	3 M Sodium formate	none	none	none

\*pH values indicated are those of the 1.0 M buffer stock solution prior to dilution with other components

