

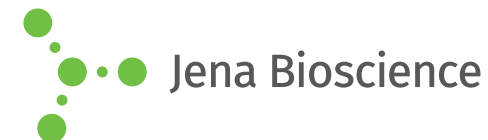


## JBScreen Classic 4

(PEG 5000 MME to 8000 based)

Cat.-No.: CS-104L

## SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
4/A1	25 % w/v Polyethylene glycol monomethyl ether 5,000	none	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
4/A2	30 % w/v Polyethylene glycol monomethyl ether 5,000	none	100 mM MES; pH 6.5	200 mM Ammonium sulfate
4/A3	3 % w/v Polyethylene glycol 6,000	none	100 mM TRIS; pH 8.5	100 mM Potassium chloride
4/A4	10 % w/v Polyethylene glycol 6,000	none	none	10 mM Magnesium chloride
4/A5	12 % w/v Polyethylene glycol 6,000	2 M Sodium chloride	none	none
4/A6	15 % w/v Polyethylene glycol 6,000	5 % w/v Glycerol	none	none
4/B1	15 % w/v Polyethylene glycol 6,000	50 mM Potassium chloride	none	10 mM Magnesium chloride
4/B2	16 % w/v Polyethylene glycol 6,000	none	none	10 mM tri-Sodium citrate
4/B3	20 % w/v Polyethylene glycol 6,000	none	50 mM Imidazole; pH 8.0	none
4/B4	25 % w/v Polyethylene glycol 6,000	none	100 mM HEPES; pH 7.5	100 mM Lithium chloride
4/B5	28 % w/v Polyethylene glycol 6,000	500 mM Lithium chloride	100 mM TRIS; pH 8.5	none
4/B6	30 % w/v Polyethylene glycol 6,000	1 M Lithium chloride	none	100 mM Sodium acetate
4/C1	33 % w/v Polyethylene glycol 6,000	none	none	10 mM tri-Sodium citrate
4/C2	2 % w/v Polyethylene glycol 8,000	500 mM Lithium sulfate	none	none
4/C3	2 % w/v Polyethylene glycol 8,000	1 M Lithium sulfate	none	none
4/C4	4 % w/v Polyethylene glycol 8,000	none	none	none
4/C5	8 % w/v Polyethylene glycol 8,000	200 mM Lithium chloride	none	50 mM Magnesium sulfate
4/C6	8 % w/v Polyethylene glycol 8,000	none	100 mM TRIS; pH 8.5	none
4/D1	10 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Zinc acetate
4/D2	10 % w/v Polyethylene glycol 8,000	none	100 mM HEPES; pH 7.5	200 mM Calcium acetate
4/D3	10 % w/v Polyethylene glycol 8,000	none	none	100 mM Sodium acetate, 50 mM Magnesium acetate
4/D4	10 % w/v Polyethylene glycol 8,000	none	none	200 mM Magnesium acetate
4/D5	10 % w/v Polyethylene glycol 8,000	10 % w/v Ethylene glycol	100 mM HEPES; pH 7.5	none
4/D6	10 % w/v Polyethylene glycol 8,000	10 % w/v Polyethylene glycol 1,000	none	none

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

