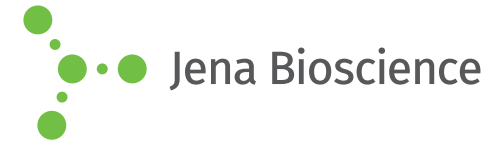




## JBScreen Classic 2

(PEG 4000 based)  
Cat.-No.: CS-102L

# SCREEN FORMULATION



Jena Bioscience

No.	Precipitant	Buffer	Additive
2/A1	4 % w/v Polyethylene glycol 4,000	100 mM Sodium acetate; pH 4.6	none
2/A2	8 % w/v Polyethylene glycol 4,000	none	none
2/A3	8 % w/v Polyethylene glycol 4,000	100 mM Sodium acetate; pH 4.6	none
2/A4	10 % w/v Polyethylene glycol 4,000	100 mM MES; pH 6.5	200 mM Magnesium chloride
2/A5	12 % w/v Polyethylene glycol 4,000	100 mM HEPES; pH 7.5	100 mM Sodium acetate
2/A6	12 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	none
2/B1	16 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
2/B2	16 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Sodium acetate
2/B3	16 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Magnesium chloride
2/B4	18 % w/v Polyethylene glycol 4,000	100 mM Sodium acetate; pH 4.6	none
2/B5	20 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
2/B6	20 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Calcium chloride
2/C1	22 % w/v Polyethylene glycol 4,000	100 mM HEPES; pH 7.5	100 mM Sodium acetate
2/C2	25 % w/v Polyethylene glycol 4,000	100 mM Sodium acetate; pH 4.6	none
2/C3	25 % w/v Polyethylene glycol 4,000	100 mM MES; pH 6.5	200 mM Magnesium chloride
2/C4	25 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Calcium chloride
2/C5	30 % w/v Polyethylene glycol 4,000	none	none
2/C6	30 % w/v Polyethylene glycol 4,000	100 mM Sodium acetate; pH 4.6	100 mM Magnesium chloride
2/D1	30 % w/v Polyethylene glycol 4,000	100 mM MES; pH 6.5	none
2/D2	30 % w/v Polyethylene glycol 4,000	100 mM HEPES; pH 7.5	200 mM Calcium chloride
2/D3	30 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
2/D4	30 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Sodium acetate
2/D5	30 % w/v Polyethylene glycol 4,000	100 mM TRIS; pH 8.5	200 mM Magnesium chloride
2/D6	35 % w/v Polyethylene glycol 4,000	none	none

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