

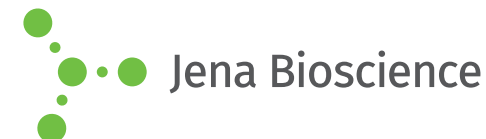


JBScreen Classic 3

(PEG 4000+ based)

Cat.-No.: CS-103L

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
3/A1	8 % w/v Polyethylene glycol 4,000	800 mM Lithium chloride	100 mM TRIS; pH 8.5	none
3/A2	10 % w/v Polyethylene glycol 4,000	20 % w/v 2-Propanol	none	none
3/A3	10 % w/v Polyethylene glycol 4,000	10 % w/v 2-Propanol	100 mM tri-Sodium citrate; pH 5.6	none
3/A4	10 % w/v Polyethylene glycol 4,000	5 % w/v 2-Propanol	100 mM HEPES; pH 7.5	none
3/A5	10 % w/v Polyethylene glycol 4,000	20 % w/v 2-Propanol	100 mM HEPES; pH 7.5	none
3/A6	12 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate
3/B1	15 % w/v Polyethylene glycol 4,000	none	none	200 mM Ammonium sulfate
3/B2	15 % w/v Polyethylene glycol 4,000	none	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium sulfate
3/B3	16 % w/v Polyethylene glycol 4,000	10 % w/v 2-Propanol	100 mM HEPES; pH 7.5	200 mM Ammonium sulfate
3/B4	20 % w/v Polyethylene glycol 4,000	none	none	200 mM Ammonium sulfate
3/B5	20 % w/v Polyethylene glycol 4,000	10 % w/v Glycerol	none	200 mM Magnesium sulfate
3/B6	20 % w/v Polyethylene glycol 4,000	5 % w/v 2-Propanol	none	100 mM tri-Sodium citrate
3/C1	20 % w/v Polyethylene glycol 4,000	20 % w/v 2-Propanol	none	100 mM tri-Sodium citrate
3/C2	20 % w/v Polyethylene glycol 4,000	none	100 mM MES; pH 6.5	600 mM Sodium chloride
3/C3	20 % w/v Polyethylene glycol 4,000	10 % w/v 2-Propanol	100 mM HEPES; pH 7.5	none
3/C4	22 % w/v Polyethylene glycol 4,000	none	none	100 mM Sodium acetate, 200 mM Ammonium sulfate
3/C5	25 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate
3/C6	25 % w/v Polyethylene glycol 4,000	none	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium sulfate
3/D1	25 % w/v Polyethylene glycol 4,000	200 mM Lithium sulfate	100 mM HEPES; pH 7.5	100 mM Sodium acetate
3/D2	25 % w/v Polyethylene glycol 4,000	8 % w/v 2-Propanol	none	100 mM Sodium acetate
3/D3	30 % w/v Polyethylene glycol 4,000	none	none	200 mM Ammonium sulfate
3/D4	30 % w/v Polyethylene glycol 4,000	none	100 mM Sodium acetate; pH 4.6	200 mM Ammonium sulfate
3/D5	30 % w/v Polyethylene glycol 4,000	none	100 mM tri-Sodium citrate; pH 5.6	100 mM Ammonium sulfate
3/D6	32 % w/v Polyethylene glycol 4,000	none	100 mM TRIS; pH 8.5	800 mM Lithium chloride

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

