



JBScreen Classic 5

based on PEG 8000 to 20000
Cat.-No.: CS-105L

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
5/A1	12 % w/v Polyethylene glycol 8,000	5 % w/v Glycerol	none	100 mM Potassium chloride
5/A2	12 % w/v Polyethylene glycol 8,000	10 % w/v Glycerol	none	500 mM Potassium chloride
5/A3	15 % w/v Polyethylene glycol 8,000	none	none	200 mM Ammonium sulfate
5/A4	15 % w/v Polyethylene glycol 8,000	500 mM Lithium sulfate	none	none
5/A5	15 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Sodium acetate
5/A6	15 % w/v Polyethylene glycol 8,000	none	none	100 mM tri-Sodium citrate, 50 mM Ammonium sulfate
5/B1	18 % w/v Polyethylene glycol 8,000	none	100 mM HEPES; pH 7.5	200 mM Calcium acetate
5/B2	18 % w/v Polyethylene glycol 8,000	2 % w/v 2-Propanol	100 mM HEPES; pH 7.5	100 mM Sodium acetate
5/B3	18 % w/v Polyethylene glycol 8,000	none	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
5/B4	20 % w/v Polyethylene glycol 8,000	none	100 mM HEPES; pH 7.5	none
5/B5	20 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Magnesium acetate
5/B6	20 % w/v Polyethylene glycol 8,000	none	100 mM CHES; pH 9.5	none
5/C1	22 % w/v Polyethylene glycol 8,000	none	100 mM MES; pH 6.5	200 mM Ammonium sulfate
5/C2	25 % w/v Polyethylene glycol 8,000	none	none	200 mM Lithium chloride
5/C3	30 % w/v Polyethylene glycol 8,000	none	none	200 mM Ammonium sulfate
5/C4	8 % w/v Polyethylene glycol 10,000	none	100 mM Sodium acetate; pH 4.6	none
5/C5	14 % w/v Polyethylene glycol 10,000	none	100 mM Imidazole; pH 8.0	none
5/C6	16 % w/v Polyethylene glycol 10,000	none	100 mM TRIS; pH 8.5	none
5/D1	18 % w/v Polyethylene glycol 10,000	20 % w/v Glycerol	100 mM TRIS; pH 8.5	100 mM Sodium chloride
5/D2	20 % w/v Polyethylene glycol 10,000	none	100 mM HEPES; pH 7.5	none
5/D3	30 % w/v Polyethylene glycol 10,000	none	100 mM TRIS; pH 8.5	none
5/D4	10 % w/v Polyethylene glycol 20,000	none	100 mM MES; pH 6.5	none
5/D5	17 % w/v Polyethylene glycol 20,000	none	100 mM TRIS; pH 8.5	100 mM Magnesium chloride





JBScreen Classic 5

based on PEG 8000 to 20000
Cat.-No.: CS-105L

SCREEN FORMULATION



5/D6	20 % w/v Polyethylene glycol 20,000	none	none	none
------	-------------------------------------	------	------	------

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