



JBScreen Membrane 2

(PEG 2000 MME to PEG 10000 based)

Cat.-No.: CS-302L

SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
C1	12 % w/v Polyethylene glycol monomethyl ether 2,000	none	50 mM TRIS; pH 8.5	500 mM Sodium chloride
C2	10 % w/v Polyethylene glycol 3,350	none	50 mM tri-Sodium citrate; pH 5.6	150 mM Sodium chloride
C3	2 % w/v Polyethylene glycol 4,000	none	50 mM TRIS; pH 7.5	none
C4	5 % w/v Polyethylene glycol 4,000	none	none	none
C5	5 % w/v Polyethylene glycol 4,000	none	none	100 mM Potassium chloride
C6	5 % w/v Polyethylene glycol 4,000	10 % w/v Glycerol	50 mM MES; pH 6.5	100 mM Sodium chloride
C7	5 % w/v Polyethylene glycol 4,000	none	50 mM Sodium Phosphate; pH 6.7	none
C8	10 % w/v Polyethylene glycol 4,000	none	50 mM TRIS; pH 8.5	500 mM Sodium chloride
C9	12 % w/v Polyethylene glycol 4,000	none	100 mM ADA; pH 6.5	100 mM Lithium sulfate
C10	12 % w/v Polyethylene glycol 4,000	none	50 mM Sodium Phosphate; pH 6.8	none
C11	12 % w/v Polyethylene glycol 4,000	20 % w/v Glycerol	50 mM MOPS; pH 7.0	500 mM Potassium chloride
C12	15 % w/v Polyethylene glycol 4,000	none	10 mM TRIS; pH 7.5	100 mM Lithium chloride
D1	20 % w/v Polyethylene glycol 4,000	none	100 mM BIS-TRIS; pH 7.0	500 mM Sodium chloride
D2	20 % w/v Polyethylene glycol 4,000	none	100 mM Sodium Phosphate; pH 7.0	500 mM Sodium chloride
D3	20 % w/v Polyethylene glycol 4,000	150 mM Zinc acetate	50 mM TRIS; pH 7.5	50 mM Zinc chloride
D4	22 % w/v Polyethylene glycol 4,000	none	50 mM TRICINE; pH 8.0	none
D5	22 % w/v Polyethylene glycol 4,000	none	50 mM TRIS; pH 8.5	500 mM Sodium chloride
D6	30 % w/v Polyethylene glycol 4,000	none	none	none
D7	10 % w/v Polyethylene glycol monomethyl ether 5,000	none	100 mM tri-Sodium citrate; pH 5.6	100 mM Magnesium acetate
D8	5 % w/v Polyethylene glycol 6,000	none	none	100 mM Magnesium sulfate
D9	10 % w/v Polyethylene glycol 6,000	150 mM Zinc acetate	50 mM TRIS; pH 7.5	50 mM Zinc chloride
D10	15 % w/v Polyethylene glycol 6,000	none	50 mM di-Sodium succinate; pH 6.5	none
D11	12 % w/v Polyethylene glycol 8,000	10 % w/v 2-Methyl-2,4-pentanediol	none	25 mM Potassium di-hydrogen phosphate
D12	8 % w/v Polyethylene glycol 10,000	none	100 mM tri-Sodium citrate; pH 5.6	100 mM Magnesium acetate

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