



## JBScreen Membrane 1

(PEG 400 to PEG 2000 MME based)

Cat.-No.: CS-301L

## SCREEN FORMULATION



No.	Precipitant	Precipitant 2	Buffer	Additive
A1	15 % w/v Polyethylene glycol 400	15 % w/v Glycerol	100 mM HEPES; pH 7.5	200 mM Calcium chloride
A2	20 % w/v Polyethylene glycol 400	100 mM Sodium chloride	100 mM tri-Sodium citrate; pH 5.6	20 mM Magnesium chloride
A3	25 % w/v Polyethylene glycol 400	none	50 mM Sodium acetate; pH 4.6	50 mM Magnesium acetate
A4	30 % w/v Polyethylene glycol 400	50 mM Sodium sulfate	50 mM TRIS; pH 8.5	50 mM Lithium sulfate
A5	48 % w/v Polyethylene glycol 400	none	100 mM HEPES; pH 7.5	200 mM Calcium chloride
A6	20 % w/v Polyethylene glycol monomethyl ether 550	none	10 mM TRIS; pH 7.5	none
A7	30 % w/v Polyethylene glycol monomethyl ether 550	none	50 mM TRIS; pH 8.5	100 mM Magnesium chloride
A8	35 % w/v Polyethylene glycol 600	none	none	none
A9	28 % w/v Polyethylene glycol 1,000	10 % w/v Glycerol	100 mM TRICINE; pH 8.0	350 mM Sodium chloride
A10	10 % w/v Polyethylene glycol 1,500	5 % w/v Ethanol	none	100 mM Magnesium chloride, 100 mM Sodium chloride
A11	30 % w/v Polyethylene glycol 1,500	none	none	none
A12	5 % w/v Polyethylene glycol 2,000	none	none	none
B1	10 % w/v Polyethylene glycol 2,000	none	100 mM TRIS; pH 8.5	500 mM Magnesium chloride
B2	15 % w/v Polyethylene glycol 2,000	none	none	none
B3	15 % w/v Polyethylene glycol 2,000	none	none	100 mM Lithium chloride
B4	15 % w/v Polyethylene glycol 2,000	none	100 mM Sodium Phosphate; pH 6.2	20 mM tri-Sodium citrate
B5	15 % w/v Polyethylene glycol 2,000	none	100 mM Sodium Phosphate; pH 6.8	500 mM Sodium chloride
B6	15 % w/v Polyethylene glycol 2,000	none	20 mM BIS-TRIS; pH 7.0	none
B7	15 % w/v Polyethylene glycol 2,000	none	50 mM HEPES; pH 7.5	100 mM Magnesium chloride
B8	20 % w/v Polyethylene glycol 2,000	2 % w/v 2-Methyl-2,4-pentanediol	100 mM TRIS; pH 8.0	300 mM Magnesium nitrate
B9	25 % w/v Polyethylene glycol 2,000	15 % w/v Glycerol	100 mM BICINE; pH 9.0	300 mM Magnesium chloride
B10	30 % w/v Polyethylene glycol 2,000	none	100 mM Sodium Phosphate; pH 6.2	250 mM Sodium chloride
B11	8 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM Sodium acetate; pH 4.6	none
B12	10 % w/v Polyethylene glycol monomethyl ether 2,000	20 % w/v Glycerol	100 mM tri-Sodium citrate; pH 5.6	3 % w/v Polyethylene glycol 200

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

