



JBScreen PACT++ 1

Cat.-No.: CS-161

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
A1	25 % w/v Polyethylene glycol 1,500	100 mM SPG buffer; pH 4.0	none
A2	25 % w/v Polyethylene glycol 1,500	100 mM SPG buffer; pH 5.0	none
A3	25 % w/v Polyethylene glycol 1,500	100 mM SPG buffer; pH 6.0	none
A4	25 % w/v Polyethylene glycol 1,500	100 mM SPG buffer; pH 7.0	none
A5	25 % w/v Polyethylene glycol 1,500	100 mM SPG buffer; pH 8.0	none
A6	25 % w/v Polyethylene glycol 1,500	100 mM SPG buffer; pH 9.0	none
A7	20 % w/v Polyethylene glycol 6,000	100 mM Sodium acetate; pH 5.0	200 mM Sodium chloride
A8	20 % w/v Polyethylene glycol 6,000	100 mM Sodium acetate; pH 5.0	200 mM Ammonium chloride
A9	20 % w/v Polyethylene glycol 6,000	100 mM Sodium acetate; pH 5.0	200 mM Lithium chloride
A10	20 % w/v Polyethylene glycol 6,000	100 mM Sodium acetate; pH 5.0	200 mM Magnesium chloride
A11	20 % w/v Polyethylene glycol 6,000	100 mM Sodium acetate; pH 5.0	200 mM Calcium chloride
A12	20 % w/v Polyethylene glycol 6,000	100 mM Sodium acetate; pH 5.0	10 mM Zinc chloride
B1	25 % w/v Polyethylene glycol 1,500	100 mM MIB buffer; pH 4.0	none
B2	25 % w/v Polyethylene glycol 1,500	100 mM MIB buffer; pH 5.0	none
B3	25 % w/v Polyethylene glycol 1,500	100 mM MIB buffer; pH 6.0	none
B4	25 % w/v Polyethylene glycol 1,500	100 mM MIB buffer; pH 7.0	none
B5	25 % w/v Polyethylene glycol 1,500	100 mM MIB buffer; pH 8.0	none
B6	25 % w/v Polyethylene glycol 1,500	100 mM MIB buffer; pH 9.0	none
B7	20 % w/v Polyethylene glycol 6,000	100 mM MES; pH 6.0	200 mM Sodium chloride
B8	20 % w/v Polyethylene glycol 6,000	100 mM MES; pH 6.0	200 mM Ammonium chloride
B9	20 % w/v Polyethylene glycol 6,000	100 mM MES; pH 6.0	200 mM Lithium chloride
B10	20 % w/v Polyethylene glycol 6,000	100 mM MES; pH 6.0	200 mM Magnesium chloride
B11	20 % w/v Polyethylene glycol 6,000	100 mM MES; pH 6.0	200 mM Calcium chloride
B12	20 % w/v Polyethylene glycol 6,000	100 mM MES; pH 6.0	10 mM Zinc chloride

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

