



JBScreen PACT++ 2

Cat.-No.: CS-162

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
C1	25 % w/v Polyethylene glycol 1,500	100 mM TBG Buffer; pH 4.0	none
C2	25 % w/v Polyethylene glycol 1,500	100 mM TBG Buffer; pH 5.0	none
C3	25 % w/v Polyethylene glycol 1,500	100 mM TBG Buffer; pH 6.0	none
C4	25 % w/v Polyethylene glycol 1,500	100 mM TBG Buffer; pH 7.0	none
C5	25 % w/v Polyethylene glycol 1,500	100 mM TBG Buffer; pH 8.0	none
C6	25 % w/v Polyethylene glycol 1,500	100 mM TBG Buffer; pH 9.0	none
C7	20 % w/v Polyethylene glycol 6,000	100 mM HEPES; pH 7.0	200 mM Sodium chloride
C8	20 % w/v Polyethylene glycol 6,000	100 mM HEPES; pH 7.0	200 mM Ammonium chloride
C9	20 % w/v Polyethylene glycol 6,000	100 mM HEPES; pH 7.0	200 mM Lithium chloride
C10	20 % w/v Polyethylene glycol 6,000	100 mM HEPES; pH 7.0	200 mM Magnesium chloride
C11	20 % w/v Polyethylene glycol 6,000	100 mM HEPES; pH 7.0	200 mM Calcium chloride
C12	20 % w/v Polyethylene glycol 6,000	100 mM HEPES; pH 7.0	10 mM Zinc chloride
D1	25 % w/v Polyethylene glycol 1,500	100 mM MMT buffer; pH 4.0	none
D2	25 % w/v Polyethylene glycol 1,500	100 mM MMT buffer; pH 5.0	none
D3	25 % w/v Polyethylene glycol 1,500	100 mM MMT buffer; pH 6.0	none
D4	25 % w/v Polyethylene glycol 1,500	100 mM MMT buffer; pH 7.0	none
D5	25 % w/v Polyethylene glycol 1,500	100 mM MMT buffer; pH 8.0	none
D6	25 % w/v Polyethylene glycol 1,500	100 mM MMT buffer; pH 9.0	none
D7	20 % w/v Polyethylene glycol 6,000	100 mM TRIS; pH 8.0	200 mM Sodium chloride
D8	20 % w/v Polyethylene glycol 6,000	100 mM TRIS; pH 8.0	200 mM Ammonium chloride
D9	20 % w/v Polyethylene glycol 6,000	100 mM TRIS; pH 8.0	200 mM Lithium chloride
D10	20 % w/v Polyethylene glycol 6,000	100 mM TRIS; pH 8.0	200 mM Magnesium chloride
D11	20 % w/v Polyethylene glycol 6,000	100 mM TRIS; pH 8.0	200 mM Calcium chloride
D12	20 % w/v Polyethylene glycol 6,000	100 mM TRIS; pH 8.0	2 mM Zinc chloride

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

