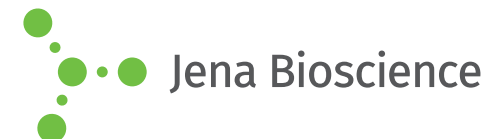




JBScreen Classic 1

based on PEG 400 to 3000
Cat.-No.: CS-101L

SCREEN FORMULATION



No.	Precipitant	Buffer	Additive
1/A1	15 % w/v Polyethylene glycol 400	100 mM Sodium acetate; pH 4.6	100 mM Calcium chloride
1/A2	15 % w/v Polyethylene glycol 400	100 mM MES; pH 6.5	none
1/A3	15 % w/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	200 mM Magnesium chloride
1/A4	15 % w/v Polyethylene glycol 400	100 mM TRIS; pH 8.5	200 mM tri-Sodium citrate
1/A5	25 % w/v Polyethylene glycol 400	100 mM Sodium acetate; pH 4.6	100 mM Magnesium chloride
1/A6	25 % w/v Polyethylene glycol 400	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
1/B1	28 % w/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	200 mM Calcium chloride
1/B2	30 % w/v Polyethylene glycol 400	100 mM Sodium acetate; pH 4.6	100 mM Calcium chloride
1/B3	30 % w/v Polyethylene glycol 400	100 mM MES; pH 6.5	100 mM Sodium acetate
1/B4	30 % w/v Polyethylene glycol 400	100 mM MES; pH 6.5	100 mM Magnesium chloride
1/B5	30 % w/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	200 mM Magnesium chloride
1/B6	30 % w/v Polyethylene glycol 400	100 mM TRIS; pH 8.5	200 mM tri-Sodium citrate
1/C1	30 % w/v Polyethylene glycol monomethyl ether 550	100 mM BICINE; pH 9.0	100 mM Sodium chloride
1/C2	25 % w/v Polyethylene glycol monomethyl ether 550	100 mM MES; pH 6.5	10 mM Zinc sulfate
1/C3	25 % w/v Polyethylene glycol 1,000	100 mM HEPES; pH 7.5	none
1/C4	30 % w/v Polyethylene glycol 1,000	100 mM TRIS; pH 8.5	none
1/C5	15 % w/v Polyethylene glycol 1,500	none	none
1/C6	20 % w/v Polyethylene glycol 1,500	100 mM HEPES; pH 7.5	none
1/D1	30 % w/v Polyethylene glycol 1,500	none	none
1/D2	20 % w/v Polyethylene glycol monomethyl ether 2,000	100 mM TRIS; pH 8.5	10 mM Nickel (II) chloride
1/D3	25 % w/v Polyethylene glycol monomethyl ether 2,000	none	none
1/D4	30 % w/v Polyethylene glycol monomethyl ether 2,000	100 mM MES; pH 6.5	100 mM Sodium acetate
1/D5	20 % w/v Polyethylene glycol 3,000	100 mM HEPES; pH 7.5	200 mM Sodium acetate
1/D6	30 % w/v Polyethylene glycol 3,000	100 mM TRIS; pH 8.5	200 mM Lithium sulfate

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

