



No.	Precipitant	Precipitant 2	Buffer	Additive
34/E1	20 % v/v Glycerol	none	none	40 mM di-Potassium hydrogen phosphate, 16 % w/v Polyethylene glycol 8,000
34/E2	15 % v/v Ethanol	none	100 mM TRIS; pH 8.0	100 mM Sodium chloride, 5 % v/v 2-Methyl-2,4-pentanediol
34/E3	40 % v/v Ethanol	none	100 mM Sodium phosphate citrate; pH 4.2	5 % w/v Polyethylene glycol 1,000
34/E4	200 mM Ammonium sulfate	none	100 mM BIS-TRIS; pH 5.5	none
34/E5	2 M Ammonium sulfate	none	100 mM Sodium acetate; pH 5.5	2 % v/v Polyethylene glycol 400
34/E6	1.6 M Ammonium sulfate	none	100 mM tri-Sodium citrate; pH 4.0	none
34/E7	2 M Lithium sulfate	none	100 mM Sodium acetate; pH 4.5	100 mM Magnesium sulfate, 5 % v/v 2-Propanol
34/E8	2 M Lithium sulfate	none	100 mM TRIS; pH 8.5	2 % v/v Polyethylene glycol 400
34/E9	2 M Lithium sulfate	none	100 mM Sodium acetate; pH 5.5	100 mM Magnesium sulfate, 5 % v/v Polyethylene glycol 400
34/E10	50 % v/v Polyethylene glycol 200	none	100 mM MES; pH 6.5	200 mM Magnesium chloride
34/E11	40 % v/v Polyethylene glycol 300	none	100 mM MES; pH 6.5	200 mM Calcium acetate
34/E12	30 % v/v Jeffamine® M-600; pH 7.0	none	100 mM HEPES; pH 7.0	none
34/F1	800 mM di-Sodium succinate; pH 7.0	none	none	none
34/F2	40 % v/v Polyethylene glycol 400	none	100 mM TRIS; pH 8.5	200 mM Lithium sulfate
34/F3	50 % v/v Polyethylene glycol 400	none	100 mM Sodium acetate; pH 4.5	200 mM Lithium sulfate
34/F4	15 % v/v Polyethylene glycol monomethyl ether 550	none	100 mM MES; pH 6.5	none
34/F5	25 % w/v Polyethylene glycol 1,500	none	100 mM SPG buffer; pH 5.5	none
34/F6	25 % w/v Polyethylene glycol 1,500	none	100 mM SPG buffer; pH 8.5	none
34/F7	25 % w/v Polyethylene glycol 1,500	none	100 mM MMT buffer; pH 6.5	none
34/F8	25 % w/v Polyethylene glycol 1,500	none	100 mM MMT buffer; pH 9.0	none
34/F9	25 % w/v Polyethylene glycol 1,500	none	100 mM MIB buffer; pH 5.0	none
34/F10	25 % w/v Polyethylene glycol 1,500	none	100 mM MIB buffer; pH 7.0	none
34/F11	12 % w/v Polyethylene glycol 1,500	none	100 mM Sodium acetate; pH 5.5	2.5 M Sodium chloride, 1.5 % v/v 2-Methyl-2,4-pentanediol
34/F12	2.4 M di-Sodium malonate	none	none	none
34/G1	30 % w/v Polyethylene glycol monomethyl ether 2,000	none	none	150 mM Potassium bromide

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



No.	Precipitant	Precipitant 2	Buffer	Additive
34/G2	10 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM Sodium acetate; pH 5.5	200 mM Ammonium sulfate
34/G3	20 % w/v Polyethylene glycol monomethyl ether 2,000	none	100 mM TRIS; pH 8.5	200 mM Trimethylamine N-oxide
34/G4	20 % w/v Polyethylene glycol 3,350	none	100 mM BIS-TRIS propane; pH 6.5	200 mM Sodium fluoride
34/G5	20 % w/v Polyethylene glycol 3,350	none	100 mM tri-Sodium citrate; pH 4.0	200 mM tri-Sodium citrate
34/G6	20 % w/v Polyethylene glycol 3,350	none	100 mM BIS-TRIS propane; pH 8.5	200 mM di-Sodium malonate
34/G7	20 % w/v Poly(acrylic acid sodium salt) 5,100	none	100 mM HEPES; pH 7.0	20 mM Magnesium chloride
34/G8	2.1 M di-Sodium DL-malate; pH 7.0	none	none	none
34/G9	800 mM di-Potassium hydrogen phosphate	800 mM Sodium di-hydrogen phosphate	100 mM HEPES; pH 7.5	200 mM Lithium chloride
34/G10	20 % w/v Polyethylene glycol 6,000	none	100 mM MES; pH 6.0	200 mM Ammonium chloride
34/G11	20 % w/v Polyethylene glycol 6,000	none	100 mM HEPES; pH 7.0	200 mM Sodium chloride
34/G12	20 % w/v Polyethylene glycol 6,000	none	100 mM TRIS; pH 8.0	200 mM Lithium chloride
34/H1	20 % w/v Polyvinylpyrrolidone K15	none	100 mM TRIS; pH 8.5	100 mM Cobalt (II) chloride
34/H2	50 % v/v Ethylene glycol	none	100 mM TRIS; pH 8.5	200 mM Magnesium chloride
34/H3	20 % w/v Polyethylene glycol 8,000	none	100 mM Imidazole; pH 6.5	3 % v/v 2-Methyl-2,4-pentanediol
34/H4	20 % w/v Polyethylene glycol 8,000	none	100 mM TRIS; pH 8.5	100 mM Magnesium chloride, 20 % v/v Polyethylene glycol 400
34/H5	20 % w/v Polyethylene glycol 8,000	none	100 mM HEPES; pH 7.5	200 mM Ammonium sulfate, 10 % v/v 2-Propanol
34/H6	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM Sodium acetate; pH 4.5	25 % w/v Polyethylene glycol 1,500
34/H7	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM Imidazole; pH 6.5	200 mM Ammonium sulfate, 10 % w/v Polyethylene glycol 3,350
34/H8	30 % v/v 2-Methyl-2,4-pentanediol	none	100 mM TRIS; pH 8.5	500 mM Sodium chloride, 8 % w/v Polyethylene glycol 8,000
34/H9	40 % v/v 2-Propanol	none	100 mM Imidazole; pH 6.5	15 % w/v Polyethylene glycol 8,000
34/H10	30 % v/v 2-Propanol	none	100 mM TRIS; pH 8.5	30 % w/v Polyethylene glycol 3,350
34/H11	17 % w/v Polyethylene glycol 10,000	none	100 mM BIS-TRIS; pH 5.5	100 mM Ammonium acetate
34/H12	15 % w/v Polyethylene glycol 20,000	none	100 mM HEPES; pH 7.0	none

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