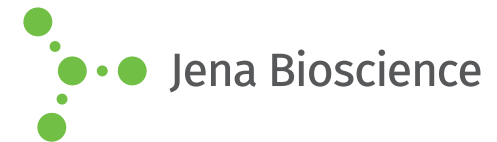




JBScreen Classic 6
(Ammonium Sulfate based)
Cat.-No.: CS-106L

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive
6/A1	500 mM Ammonium sulfate	1 M Lithium sulfate	none	100 mM tri-Sodium citrate
6/A2	1 M Ammonium sulfate	none	none	none
6/A3	1 M Ammonium sulfate	none	100 mM Sodium acetate; pH 4.6	none
6/A4	1 M Ammonium sulfate	2 % w/v Polyethylene glycol 400	100 mM HEPES; pH 7.5	none
6/A5	1 M Ammonium sulfate	none	100 mM TRIS; pH 8.5	none
6/A6	1.2 M Ammonium sulfate	3 % w/v 2-Propanol	none	50 mM tri-Sodium citrate
6/B1	1.5 M Ammonium sulfate	15 % w/v Glycerol	100 mM TRIS; pH 8.5	none
6/B2	1.6 M Ammonium sulfate	500 mM Lithium chloride	none	none
6/B3	1.6 M Ammonium sulfate	1 M Lithium sulfate	none	none
6/B4	1.6 M Ammonium sulfate	none	100 mM HEPES; pH 7.5	200 mM Sodium chloride
6/B5	1.6 M Ammonium sulfate	2 % w/v Polyethylene glycol 1,000	100 mM HEPES; pH 7.5	none
6/B6	1.8 M Ammonium sulfate	none	100 mM MES; pH 6.5	none
6/C1	2 M Ammonium sulfate	2 M Sodium chloride	none	none
6/C2	2 M Ammonium sulfate	none	100 mM Sodium acetate; pH 4.6	none
6/C3	2 M Ammonium sulfate	5 % w/v Polyethylene glycol 400	100 mM MES; pH 6.5	none
6/C4	2 M Ammonium sulfate	none	100 mM TRIS; pH 8.5	none
6/C5	2.2 M Ammonium sulfate	none	none	none
6/C6	2.2 M Ammonium sulfate	20 % w/v Glycerol	none	none
6/D1	2.4 M Ammonium sulfate	none	none	100 mM tri-Sodium citrate
6/D2	3 M Ammonium sulfate	1 % w/v 2-Methyl-2,4-pentanediol	none	none
6/D3	3 M Ammonium sulfate	10 % w/v Glycerol	none	none
6/D4	3.5 M Ammonium sulfate	none	100 mM HEPES; pH 7.5	none
6/D5	3.5 M Ammonium sulfate	1 % w/v 2-Methyl-2,4-pentanediol	100 mM MES; pH 6.5	none
6/D6	3.5 M Ammonium sulfate	none	none	none

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