



XP Screen

Crystallization Screen for improved Crystal Quality and Phasing

Cat.-No.: CS-350

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive 1	Additive 2
A1	25.00 % v/v Ethylene glycol	None	None		1 mM TEW
A2	12.00 % v/v Glycerol	1.5 M Ammonium sulfate	100 mM TRIS; pH 8.5		1 mM TEW
A3	100 mM 1,6-Hexanediol	None	100 mM Sodium acetate; pH 4.6	5.00 mM Cobalt (II) chloride	1 mM TEW
A4	2.5 M 1,6-Hexanediol	None	100 mM tri-Sodium citrate; pH 5.6		1 mM TEW
A5	1.00 M 1,6-Hexanediol	None	100 mM HEPES; pH 7.5	10.00 mM Magnesium chloride	1 mM TEW
A6	30.00 % v/v 2-Methyl-2,4-pentanediol	None	100 mM Sodium acetate; pH 4.6	10.00 mM Sodium chloride	1 mM TEW
A7	30.00 % v/v 2-Methyl-2,4-pentanediol	None	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium acetate	1 mM TEW
A8	10.00 % v/v 2-Methyl-2,4-pentanediol	None	100 mM Sodium acetate; pH 4.6	5.00 mM Calcium chloride	1 mM TEW
A9	30.00 % v/v 2-Methyl-2,4-pentanediol	500 mM Ammonium sulfate	100 mM HEPES; pH 7.5		1 mM TEW
A10	30.00 % v/v 2-Methyl-2,4-pentanediol	None	100 mM HEPES; pH 7.5	200 mM tri-Sodium citrate	1 mM TEW
A11	50.00 % v/v 2-Methyl-2,4-pentanediol	None	100 mM TRIS; pH 8.5	200 mM Ammonium di-hydrogen phosphate	1 mM TEW
A12	50.00 % v/v 2-Methyl-2,4-pentanediol	None	100 mM HEPES; pH 7.5		1 mM TEW
B1	2.00 % v/v Polyethylene glycol 400	None	100 mM tri-Sodium citrate; pH 5.6	500 mM Sodium chloride	1 mM TEW
B2	2.00 % v/v Polyethylene glycol 400	2.00 M Ammonium sulfate	100 mM HEPES; pH 7.5		1 mM TEW
B3	10.00 % v/v Polyethylene glycol 400	None	100 mM HEPES; pH 7.5	10.00 mM Calcium chloride	1 mM TEW
B4	30.00 % v/v Polyethylene glycol 400	None	100 mM tri-Sodium citrate; pH 5.6		1 mM TEW
B5	30.00 % v/v Polyethylene glycol 400	None	100 mM HEPES; pH 7.5	200 mM Magnesium chloride	1 mM TEW
B6	10.00 % v/v Polyethylene glycol 400	None	100 mM Sodium acetate; pH 4.6	10.00 mM Calcium chloride	1 mM TEW
B7	20.00 % v/v Polyethylene glycol monomethyl ether 550	None	100 mM HEPES; pH 7.5	100 mM Sodium chloride	1 mM TEW
B8	25.00 % v/v Polyethylene glycol monomethyl ether 550	None	100 mM MES; pH 6.5		1 mM TEW
B9	10.00 % w/v Polyethylene glycol 1,000	10.00 % w/v Polyethylene glycol 8,000	100 mM tri-Sodium citrate; pH 5.6		1 mM TEW
B10	30.00 % w/v Polyethylene glycol 1,500	None	100 mM HEPES; pH 7.5		1 mM TEW
B11	5.00 % w/v Polyethylene glycol monomethyl ether 2,000	None	100 mM TRIS; pH 8.5		1 mM TEW
B12	20.00 % w/v Polyethylene glycol monomethyl ether 2,000	None	None		1 mM TEW

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



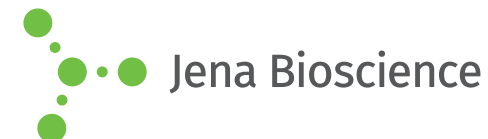


XP Screen

Crystallization Screen for improved Crystal Quality and Phasing

Cat.-No.: CS-350

SCREEN FORMULATION



Jena Bioscience

No.	Precipitant 1	Precipitant 2	Buffer	Additive 1	Additive 2
C1	30.00 % w/v Polyethylene glycol monomethyl ether 2,000	None	100 mM Sodium acetate; pH 4.6	10.00 mM Ammonium sulfate	1 mM TEW
C2	8.00 % w/v Polyethylene glycol 4,000	None	100 mM Sodium acetate; pH 4.6		1 mM TEW
C3	10.00 % w/v Polyethylene glycol 4,000	5.00 % v/v 2-Propanol	100 mM tri-Sodium citrate; pH 5.6		1 mM TEW
C4	20.00 % w/v Polyethylene glycol 4,000	10.00 % v/v 2-Propanol	100 mM HEPES; pH 7.5		1 mM TEW
C5	10.00 % w/v Polyethylene glycol 4,000	None	100 mM Sodium acetate; pH 4.6	10.00 mM Ammonium sulfate	1 mM TEW
C6	30.00 % w/v Polyethylene glycol 4,000	None	None	200 mM Ammonium sulfate	1 mM TEW
C7	15.00 % w/v Polyethylene glycol 4,000	None	100 mM Sodium acetate; pH 4.6	10.00 mM Ammonium acetate	1 mM TEW
C8	30.00 % w/v Polyethylene glycol 4,000	None	100 mM tri-Sodium citrate; pH 5.6	200 mM Ammonium acetate	1 mM TEW
C9	30.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	200 mM Sodium acetate	1 mM TEW
C10	30.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	200 mM Lithium sulfate	1 mM TEW
C11	30.00 % w/v Polyethylene glycol 4,000	None	100 mM TRIS; pH 8.5	10.00 mM Magnesium chloride	1 mM TEW
C12	10.00 % w/v Polyethylene glycol monomethyl ether 5,000	None	100 mM MES; pH 6.5	10.00 mM Ammonium sulfate	1 mM TEW
D1	10.00 % w/v Polyethylene glycol 6,000	None	100 mM tri-Sodium citrate; pH 5.6	100 mM Sodium chloride	1 mM TEW
D2	10.00 % w/v Polyethylene glycol 6,000	5.00 % v/v 2-Methyl-2,4-pentanediol	100 mM HEPES; pH 7.5		1 mM TEW
D3	8.00 % w/v Polyethylene glycol 8,000	None	100 mM TRIS; pH 8.5		1 mM TEW
D4	10.00 % w/v Polyethylene glycol 8,000	8.00 % v/v Ethylene glycol	100 mM HEPES; pH 7.5		1 mM TEW
D5	15.00 % w/v Polyethylene glycol 8,000	500 mM Lithium sulfate	100 mM Sodium acetate; pH 4.6		1 mM TEW
D6	10.00 % w/v Polyethylene glycol 8,000	None	100 mM MES; pH 6.5	10.00 mM Calcium acetate	1 mM TEW
D7	18.00 % w/v Polyethylene glycol 8,000	None	100 mM MES; pH 6.5		1 mM TEW
D8	10.00 % w/v Polyethylene glycol 8,000	None	None	10.00 mM Potassium di-hydrogen phosphate	1 mM TEW
D9	20.00 % w/v Polyethylene glycol 8,000	None	100 mM MES; pH 6.5	200 mM Magnesium acetate	1 mM TEW
D10	10.00 % w/v Polyethylene glycol 8,000	None	100 mM MES; pH 6.5	5.00 mM Sodium acetate	1 mM TEW
D11	30.00 % w/v Polyethylene glycol 8,000	None	None	200 mM Ammonium sulfate	1 mM TEW
D12	10.00 % w/v Polyethylene glycol 8,000	None	100 mM MES; pH 6.5	5.00 mM Ammonium sulfate	1 mM TEW

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



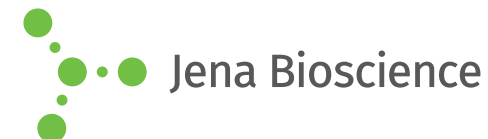


XP Screen

Crystallization Screen for improved Crystal Quality and Phasing

Cat.-No.: CS-350

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive 1	Additive 2
E1	10.00 % w/v Polyethylene glycol 10,000	2.00 % v/v 1,4-Dioxane	100 mM tri-Sodium citrate; pH 5.6		1 mM TEW
E2	20.00 % w/v Polyethylene glycol 10,000	None	100 mM HEPES; pH 7.5		1 mM TEW
E3	12.00 % w/v Polyethylene glycol 20,000	None	100 mM MES; pH 6.5		1 mM TEW
E4	5.00 % v/v 2-Propanol	2.00 M Ammonium sulfate	None		1 mM TEW
E5	20.00 % v/v 2-Propanol	None	100 mM HEPES; pH 7.5	200 mM tri-Sodium citrate	1 mM TEW
E6	20.00 % v/v 2-Propanol	None	100 mM Sodium acetate; pH 4.6		1 mM TEW
E7	30.00 % v/v 2-Propanol	None	100 mM HEPES; pH 7.5	200 mM Magnesium chloride	1 mM TEW
E8	10.00 % v/v 2-Propanol	None	100 mM TRIS; pH 8.5	10.00 mM Ammonium acetate	1 mM TEW
E9	10.00 % v/v 1,4-Dioxane	1.6 M Ammonium sulfate	100 mM MES; pH 6.5		1 mM TEW
E10	35.00 % v/v 1,4-Dioxane	None	None		1 mM TEW
E11	10.00 % v/v Ethanol	None	None	100 mM Sodium chloride	1 mM TEW
E12	20.00 % v/v Ethanol	None	100 mM TRIS; pH 8.5		1 mM TEW
F1	25.00 % v/v 2-Methyl-2-propanol	None	100 mM TRIS; pH 8.5		1 mM TEW
F2	10.00 % v/v 2-Methyl-2-propanol	None	100 mM HEPES; pH 7.5		1 mM TEW
F3	1.00 M Imidazole; pH 7.0	None	None		1 mM TEW
F4	100 mM Lithium sulfate	None	100 mM TRIS; pH 8.5		1 mM TEW
F5	1.5 M Lithium sulfate	None	100 mM HEPES; pH 7.5		1 mM TEW
F6	400 mM Potassium Sodium tartrate	None	None		1 mM TEW
F7	400 mM Potassium Sodium tartrate	None	100 mM HEPES; pH 7.5		1 mM TEW
F8	1.4 M tri-Sodium citrate	None	100 mM HEPES; pH 7.5		1 mM TEW
F9	1.6 M tri-Sodium citrate; pH 6.5	None	None		1 mM TEW
F10	10.00 % v/v Jeffamine® M-600; pH 7.0	None	100 mM tri-Sodium citrate; pH 5.6	10.00 mM Iron (III) chloride	1 mM TEW
F11	20.00 % v/v Jeffamine® M-600; pH 7.0	None	100 mM HEPES; pH 7.5		1 mM TEW
F12	20.00 % v/v Jeffamine® M-600; pH 7.0	None	100 mM MES; pH 6.5	10.00 mM Cesium chloride	1 mM TEW

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





XP Screen

Crystallization Screen for improved Crystal Quality and Phasing

Cat.-No.: CS-350

SCREEN FORMULATION



No.	Precipitant 1	Precipitant 2	Buffer	Additive 1	Additive 2
G1	100 mM Potassium di-hydrogen phosphate	100 mM Sodium di-hydrogen phosphate	100 mM tri-Sodium citrate; pH 5.6		1 mM TEW
G2	400 mM Ammonium di-hydrogen phosphate	None	None		1 mM TEW
G3	1.00 M Ammonium di-hydrogen phosphate	None	100 mM tri-Sodium citrate; pH 5.6		1 mM TEW
G4	2.00 M Ammonium di-hydrogen phosphate	None	100 mM TRIS; pH 8.5		1 mM TEW
G5	2.00 M Ammonium formate	None	100 mM Sodium acetate; pH 4.6		1 mM TEW
G6	4.00 M Ammonium formate	None	100 mM HEPES; pH 7.5		1 mM TEW
G7	2.00 M Ammonium formate	None	None		1 mM TEW
G8	500 mM Ammonium sulfate	1.00 M Lithium sulfate	100 mM tri-Sodium citrate; pH 5.6		1 mM TEW
G9	1.6 M Ammonium sulfate	None	100 mM HEPES; pH 7.5	100 mM Sodium chloride	1 mM TEW
G10	1.8 M Ammonium sulfate	None	100 mM MES; pH 6.5	10.00 mM Cobalt (II) chloride	1 mM TEW
G11	2.00 M Ammonium sulfate	None	100 mM TRIS; pH 8.5		1 mM TEW
G12	2.00 M Ammonium sulfate	None	None		1 mM TEW
H1	2.00 M Ammonium sulfate	None	100 mM Sodium acetate; pH 4.6		1 mM TEW
H2	2.00 M Ammonium sulfate	None	100 mM tri-Sodium citrate; pH 5.6	200 mM Potassium Sodium tartrate	1 mM TEW
H3	200 mM Magnesium formate	None	None		1 mM TEW
H4	1.6 M Magnesium sulfate	None	100 mM MES; pH 6.5		1 mM TEW
H5	2.00 M Magnesium chloride	None	100 mM BICINE; pH 9.0		1 mM TEW
H6	1.00 M Sodium acetate	None	100 mM Imidazole; pH 6.5		1 mM TEW
H7	100 mM Sodium acetate	None	100 mM tri-Sodium citrate; pH 5.6	5.00 mM Cadmium sulfate	1 mM TEW
H8	1.4 M Sodium acetate	None	100 mM MES; pH 6.5		1 mM TEW
H9	100 mM Sodium chloride	None	None	10.00 mM Magnesium chloride	1 mM TEW
H10	2.00 M Sodium chloride	None	100 mM Sodium acetate; pH 4.6		1 mM TEW
H11	2.00 M Sodium chloride	None	100 mM MES; pH 6.5	100 mM Sodium di-hydrogen phosphate, 100 mM Potassium di-hydrogen phosphate	1 mM TEW
H12	4.3 M Sodium chloride	None	100 mM HEPES; pH 7.5		1 mM TEW

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

