



Cat# EBS-BLCP

Cubic™ lipid-compatible crystallization screen - technical sheet

Formulations: (Patent No. 6,821,487)

	<u>crystallant</u>	<u>buffer (0.1 M)</u>	<u>salt (0.2 M)</u>
1	10% (v/v) 2-propanol	Tris pH 7.0	Zn(OAc) ₂
2	10% (v/v) 2-propanol	acetate pH 4.5	Ca(OAc) ₂
3	10% (v/v) 2-propanol	citrate pH 5.5	Li ₂ SO ₄
4	10% (v/v) 2-propanol	Tris pH 7.0	Ca(OAc) ₂
5	10% (v/v) 2-propanol	acetate pH 4.5	NaCl
6	10% (v/v) 2-propanol	citrate pH 5.5	MgCl ₂
7	10% (w/v) PEG-8000	HEPES pH 7.5	NaCl
8	10% (w/v) PEG-8000	MES pH 6.0	Zn(OAc) ₂
9	10% (w/v) PEG-8000	Na/K phosphate pH 6.2	Li ₂ SO ₄
10	10% (w/v) PEG-8000	acetate pH 4.5	Ca(OAc) ₂
11	10% (w/v) PEG-8000	Tris pH 8.5	MgCl ₂
12	10% (w/v) PEG-8000	cacodylate pH 6.5	NaCl
13	10% (w/v) PEG-8000	citrate pH 5.5	Li ₂ SO ₄
14	10% (w/v) PEG-8000	HEPES pH 7.5	MgCl ₂
15	20% (w/v) PEG-2000 MME	HEPES pH 7.5	NaCl
16	20% (w/v) PEG-2000 MME	MES pH 6.0	Zn(OAc) ₂
17	20% (w/v) PEG-2000 MME	Na/K phosphate pH 6.2	Li ₂ SO ₄
18	20% (w/v) PEG-2000 MME	acetate pH 4.5	Ca(OAc) ₂
19	20% (w/v) PEG-2000 MME	Tris pH 8.5	MgCl ₂
20	20% (w/v) PEG-2000 MME	cacodylate pH 6.5	Li ₂ SO ₄
21	20% (w/v) PEG-2000 MME	HEPES pH 7.5	Li ₂ SO ₄
22	20% (v/v) 1,4-butanediol	Tris pH 8.5	Li ₂ SO ₄
23	20% (v/v) 1,4-butanediol	cacodylate pH 6.5	Ca(OAc) ₂
24	20% (v/v) 1,4-butanediol	citrate pH 5.5	MgCl ₂
25	20% (v/v) 1,4-butanediol	Tris pH 7.0	Zn(OAc) ₂
26	20% (v/v) 1,4-butanediol	Tris pH 8.5	MgCl ₂
27	20% (v/v) 1,4-butanediol	citrate pH 5.5	NaCl
28	20% (w/v) PEG-1000	HEPES pH 7.5	NaCl
29	20% (w/v) PEG-1000	MES pH 6.0	Zn(OAc) ₂
30	20% (w/v) PEG-1000	Tris pH 7.0	Li ₂ SO ₄
31	20% (w/v) PEG-1000	HEPES pH 7.5	MgCl ₂
32	20% (w/v) PEG-1000	MES pH 6.0	NaCl
33	2.5 M NaCl	HEPES pH 7.5	
34	2.5 M NaCl	MES pH 6.0	Zn(OAc) ₂
35	2.5 M NaCl	Tris pH 8.5	Li ₂ SO ₄
36	2.5 M NaCl	cacodylate pH 6.5	Ca(OAc) ₂
37	2.5 M NaCl	citrate pH 5.5	MgCl ₂
38	2.5 M NaCl	cacodylate pH 6.5	Li ₂ SO ₄
39	30% (w/v) PEG-400	Tris pH 7.0	Zn(OAc) ₂
40	30% (w/v) PEG-400	Imidazole pH 8.0	Li ₂ SO ₄
41	30% (w/v) PEG-400	citrate pH 5.5	MgCl ₂
42	30% (w/v) PEG-400	citrate pH 5.5	Li ₂ SO ₄
43	30% (w/v) PEG-400	Imidazole pH 8.0	Zn(OAc) ₂
44	30% (w/v) PEG-400	Tris pH 7.0	NaCl

45	15% (v/v) Ethanol	Na/K phosphate pH 6.2	NaCl
46	15% (v/v) Ethanol	acetate pH 4.5	Zn(OAc) ₂
47	15% (v/v) Ethanol	cacodylate pH 6.5	Ca(OAc) ₂
48	15% (v/v) Ethanol	acetate pH 4.5	MgCl ₂

49	15% (v/v) Ethanol	cacodylate pH 6.5	NaCl
50	10% (w/v) PEG-3000	HEPES pH 7.5	NaCl
51	10% (w/v) PEG-3000	MES pH 6.0	Zn(OAc) ₂
52	10% (w/v) PEG-3000	Tris pH 8.5	Ca(OAc) ₂
53	10% (w/v) PEG-3000	citrate pH 5.5	MgCl ₂
54	10% (w/v) PEG-3000	Tris pH 8.5	Li ₂ SO ₄
55	10% (w/v) PEG-3000	MES pH 6.0	Li ₂ SO ₄
56	1.0 M (NH ₄) ₂ HPO ₄	HEPES pH 7.5	NaCl
57	1.0 M (NH ₄) ₂ HPO ₄	cacodylate pH 6.5	NaCl
58	1.26 M (NH ₄) ₂ SO ₄	Na/K phosphate pH 6.2	Li ₂ SO ₄
59	1.26 M (NH ₄) ₂ SO ₄	Tris pH 7.0	MgCl ₂
60	1.26 M (NH ₄) ₂ SO ₄	Imidazole pH 8.0	Li ₂ SO ₄
61	1.26 M (NH ₄) ₂ SO ₄	citrate pH 5.5	NaCl
62	20% (w/v) PEG-8000	HEPES pH 7.5	NaCl
63	20% (w/v) PEG-8000	Tris pH 7.0	Zn(OAc) ₂
64	20% (w/v) PEG-8000	Imidazole pH 8.0	Li ₂ SO ₄
65	20% (w/v) PEG-8000	acetate pH 4.5	MgCl ₂
66	20% (w/v) PEG-8000	cacodylate pH 6.5	NaCl
67	20% (w/v) PEG-8000	Tris pH 7.0	MgCl ₂
68	1.0 M Sodium citrate	HEPES pH 7.5	NaCl
69	1.0 M Sodium citrate	MES pH 6.0	Zn(OAc) ₂
70	1.0 M Sodium citrate	Imidazole pH 8.0	Li ₂ SO ₄
71	1.0 M Sodium citrate	acetate pH 4.5	MgCl ₂
72	1.0 M Sodium citrate	Tris pH 8.5	Li ₂ SO ₄
73	1.0 M Sodium citrate	MES pH 6.0	MgCl ₂
74	10% (v/v) 2-propanol	acetate pH 4.5	Li ₂ SO ₄
75	10% (v/v) 2-propanol	citrate pH 5.5	NaCl
76	10% (w/v) PEG-8000	acetate pH 4.5	MgCl ₂
77	10% (w/v) PEG-8000	Tris pH 8.5	NaCl
78	20% (w/v) PEG-2000 MME	acetate pH 4.5	Zn(OAc) ₂
79	20% (w/v) PEG-2000 MME	Tris pH 8.5	Li ₂ SO ₄
80	20% (v/v) 1,4-butanediol	Tris pH 8.5	NaCl
81	20% (v/v) 1,4-butanediol	cacodylate pH 6.5	Li ₂ SO ₄
82	20% (w/v) PEG-1000	HEPES pH 7.5	Li ₂ SO ₄
83	20% (w/v) PEG-1000	MES pH 6.0	MgCl ₂
84	2.5 M NaCl	MES pH 6.0	Li ₂ SO ₄
85	30% (w/v) PEG-8000	Tris pH 8.5	Li ₂ SO ₄
86	30% (w/v) PEG-8000	cacodylate pH 6.5	Ca(OAc) ₂
87	30% (w/v) PEG-400	MES pH 6.0	MgCl ₂
88	30% (w/v) PEG-400	Tris pH 7.0	Li ₂ SO ₄
89	10% (w/v) PEG-3000	HEPES pH 7.5	Li ₂ SO ₄
90	10% (w/v) PEG-3000	MES pH 6.0	MgCl ₂
91	1.0 M Sodium citrate	Imidazole pH 8.0	MgCl ₂
92	1.0 M Sodium citrate	Tris pH 8.5	NaCl
93	2.5M Sodium Malonate	Tris pH 7.0	
94	2.5M Sodium Malonate	acetate pH 4.5	
95	2.5M Sodium Malonate	Tris pH 8.5	
96	2.5M Sodium Malonate	Imidazole pH 8.0	

All formulations are made with ultrapure ASTM Type I water and sterile-filtered stock solutions. Store at 4-25 °C.

© 2004 Emerald BioSystems, Inc., (888) 780-8535, FAX (206) 780-8549, www.emeraldbiosystems.com