

Cryo I sparse matrix crystallization screen - technical sheet

Formulations: (Patent No. 6,267,935)

	<u>crystallant</u>	<u>buffer (0.1 M)</u>	<u>additive(s)</u>	
1	40% (v/v) 2-methyl-2,4-pentanediol	phosphate-citrate pH 4.2	none	1
2	40% (v/v) ethylene glycol	acetate pH 4.5	none	2
3	50% (v/v) PEG-200	citrate pH 5.5	none	3
4	40% (v/v) PEG-300	HEPES pH 7.5	0.2 M NaCl	4
5	40% (v/v) PEG-400	citrate pH 5.5	0.2 M MgCl ₂	5
6	40% (v/v) PEG-600	cacodylate pH 6.5	0.2 M Ca(OAc) ₂	6
7	40% (v/v) ethanol	Tris pH 8.5	0.05 M MgCl ₂	7
8	35% (v/v) 2-ethoxyethanol	cacodylate pH 6.5	none	8
9	35% (v/v) 2-propanol	phosphate-citrate pH 4.2	none	9
10	45% (v/v) glycerol	imidazole pH 8.0	none	10
11	35% (v/v) 2-methyl-2,4-pentanediol	Tris pH 8.5	0.2 M (NH ₄) ₂ SO ₄	11
12	50% (v/v) ethylene glycol	acetate pH 4.5	5% (w/v) PEG-1000	12
13	30% (v/v) PEG-200	MES pH 6.0	5% (w/v) PEG-3000	13
14	20% (v/v) PEG-300	phosphate-citrate pH 4.2	0.2 M (NH ₄) ₂ SO ₄ , 10% (v/v) glycerol	14
15	50% (v/v) PEG-400	CHES pH 9.5	0.2 M NaCl	15
16	30% (v/v) PEG-600	MES pH 6.0	5% (w/v) PEG-1000, 10% (v/v) glycerol	16
17	40% (v/v) 1, 2-propanediol	HEPES pH 7.5	none	17
18	35% (v/v) 2-ethoxyethanol	imidazole pH 8.0	0.05 M Ca(OAc) ₂	18
19	35% (v/v) 2-propanol	Tris pH 8.5	none	19
20	30% (v/v) 1,2-propanediol	citrate pH 5.5	20% (v/v) 2-methyl-2,4-pentanediol	20
21	40% (v/v) 1,2-propanediol	acetate pH 4.5	0.05 M Ca(OAc) ₂	21
22	40% (v/v) ethylene glycol	Na/K phosphate pH 6.2	none	22
23	40% (v/v) 2-methyl-2,4-pentanediol	Tris pH 7.0	0.2 M (NH ₄) ₂ SO ₄	23
24	40% (v/v) PEG-400	Na/K phosphate pH 6.2	0.2 M NaCl	24
25	30% (v/v) PEG-200	Tris pH 8.5	0.2 M (NH ₄) ₂ HPO ₄	25
26	40% (v/v) PEG-300	CHES pH 9.5	0.2 M NaCl	26
27	30% (v/v) PEG-400	CAPS pH 10.5	0.5 M (NH ₄) ₂ SO ₄ , 10% (v/v) glycerol	27
28	30% (v/v) PEG-600	HEPES pH 7.5	0.05 M Li ₂ SO ₄ , 10% (v/v) glycerol	28
29	40% (v/v) PEG-300	CHES pH 9.5	0.2 M sodium citrate	29
30	35% (v/v) 2-ethoxyethanol	citrate pH 5.5	none	30
31	35% (v/v) 2-propanol	citrate pH 5.5	5% (w/v) PEG-1000	31
32	40% (v/v) 1,2-propanediol	CHES pH 9.5	0.2 M sodium citrate	32
33	25% (v/v) 1, 2-propanediol	imidazole pH 8.0	0.2 M Zn(OAc) ₂ , 10% (v/v) glycerol	33
34	40% (v/v) 2-methyl-2,4-pentanediol	imidazole pH 8.0	0.2 M MgCl ₂	34
35	40% (v/v) ethylene glycol	HEPES pH 7.5	5% (w/v) PEG-3000	35
36	50% (v/v) PEG-200	Tris pH 7.0	0.05 M Li ₂ SO ₄	36
37	40% (v/v) PEG-300	cacodylate pH 6.5	0.2 M Ca(OAc) ₂	37
38	40% (v/v) PEG-400	Tris pH 8.5	0.2 M Li ₂ SO ₄	38
39	40% (v/v) PEG-600	phosphate-citrate pH 4.2	none	39
40	40% (v/v) ethanol	phosphate-citrate pH 4.2	5% (w/v) PEG-1000	40
41	25% (v/v) 1, 2-propanediol	phosphate-citrate pH 4.2	5% (w/v) PEG-3000, 10% (v/v) glycerol	41
42	40% (v/v) ethylene glycol	Tris pH 7.0	none	42
43	50% (v/v) ethylene glycol	Tris pH 8.5	0.2 M MgCl ₂	43
44	50% (v/v) PEG-200	cacodylate pH 6.5	0.2 M Zn(OAc) ₂	44
45	20% (v/v) PEG-300	Tris pH 8.5	5% (w/v) PEG-8000, 10% (v/v) glycerol	45
46	40% (v/v) PEG-400	MES pH 6.0	5% (w/v) PEG-3000	46
47	50% (v/v) PEG-400	acetate pH 4.5	0.2 M Li ₂ SO ₄	47
48	40% (v/v) PEG-600	imidazole pH 8.0	0.2 M Zn(OAc) ₂	48

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

Cryo II sparse matrix crystallization screen - technical sheet

Formulations: (Patent No. 6,267,935)

	<u>crystallant</u>	<u>buffer (0.1 M)</u>	<u>additive(s)</u>	
1	40% (v/v) 2-methyl-2,4-pentanediol	cacodylate pH 6.5	5% (w/v) PEG-8000	1
2	50% (v/v) PEG-200	CHES pH 9.5	none	2
3	40% (v/v) ethylene glycol	phosphate-citrate pH 4.2	0.2 M (NH ₄) ₂ SO ₄	3
4	40% (v/v) PEG-400	HEPES pH 7.5	0.2 M Ca(OAc) ₂	4
5	40% (v/v) PEG-300	Tris pH 7.0	5% (w/v) PEG-1000	5
6	30% (v/v) PEG-600	cacodylate pH 6.5	1 M NaCl, 10% (v/v) glycerol	6
7	40% (v/v) ethanol	Tris pH 7.0	none	7
8	35% (v/v) 2-ethoxyethanol	Na/K phosphate pH 6.2	0.2 M NaCl	8
9	35% (v/v) 2-propanol	imidazole pH 8.0	0.05 M Zn(OAc) ₂	9
10	40% (v/v) 1,2-propanediol	acetate pH 4.5	none	10
11	25% (v/v) 1, 2-propanediol	Na/K phosphate pH 6.2	10% (v/v) glycerol	11
12	40% (v/v) 1,2-propanediol	citrate pH 5.5	0.2 M NaCl	12
13	35% (v/v) 2-methyl-2,4-pentanediol	cacodylate pH 6.5	0.05 M Zn(OAc) ₂	13
14	40% (v/v) ethylene glycol	imidazole pH 8.0	0.2 M Ca(OAc) ₂	14
15	50% (v/v) PEG-200	Na/K phosphate pH 6.2	0.2 M NaCl	15
16	20% (v/v) PEG-300	imidazole pH 8.0	1 M (NH ₄) ₂ SO ₄ , 10% (v/v) glycerol	16
17	50% (v/v) PEG-400	MES pH 6.0	none	17
18	40% (v/v) PEG-300	phosphate-citrate pH 4.2	none	18
19	40% (v/v) PEG-600	acetate pH 4.5	0.2 M MgCl ₂	19
20	50% (v/v) ethylene glycol	CHES pH 9.5	0.5 M K/Na tartrate	20
21	35% (v/v) 2-ethoxyethanol	Tris pH 8.5	0.2 M Li ₂ SO ₄	21
22	35% (v/v) 2-propanol	cacodylate pH 6.5	0.2 M MgCl ₂	22
23	30% (v/v) 1,2-propanediol	HEPES pH 7.5	20% (v/v) PEG-400	23
24	25% (v/v) 1, 2-propanediol	Tris pH 8.5	0.2 M MgCl ₂ , 10% (v/v) glycerol	24
25	40% (v/v) 2-methyl-2,4-pentanediol	CAPS pH 10.5	none	25
26	40% (v/v) ethylene glycol	MES pH 6.0	0.2 M Zn(OAc) ₂	26
27	50% (v/v) PEG-200	Tris pH 7.0	none	27
28	40% (v/v) PEG-300	imidazole pH 8.0	0.2 M Zn(OAc) ₂	28
29	30% (v/v) PEG-400	HEPES pH 7.5	5% (w/v) PEG-3000, 10% (v/v) glycerol	29
30	40% (v/v) PEG-600	citrate pH 5.5	none	30
31	40% (v/v) PEG-600	CHES pH 9.5	none	31
32	35% (v/v) 2-propanol	acetate pH 4.5	none	32
33	45% (v/v) glycerol	cacodylate pH 6.5	0.2 M Ca(OAc) ₂	33
34	25% (v/v) 1, 2-propanediol	Tris pH 7.0	0.2 M (NH ₄) ₂ SO ₄ , 10% (v/v) glycerol	34
35	40% (v/v) 2-methyl-2,4-pentanediol	citrate pH 5.5	none	35
36	50% (v/v) PEG-200	cacodylate pH 6.5	0.2 M MgCl ₂	36
37	50% (v/v) ethylene glycol	imidazole pH 8.0	none	37
38	40% (v/v) PEG-400	acetate pH 4.5	none	38
39	30% (v/v) PEG-600	Tris pH 7.0	0.5 M (NH ₄) ₂ SO ₄ , 10% (v/v) glycerol	39
40	40% (v/v) 2-methyl-2,4-pentanediol	CHES pH 9.5	none	40
41	50% (v/v) ethylene glycol	HEPES pH 7.5	0.2 M Li ₂ SO ₄	41
42	30% (v/v) PEG-200	acetate pH 4.5	0.1 M NaCl	42
43	40% (v/v) PEG-400	imidazole pH 8.0	none	43
44	35% (v/v) 2-methyl-2,4-pentanediol	acetate pH 4.5	10% (v/v) glycerol	44
45	40% (v/v) PEG-300	acetate pH 4.5	0.2 M NaCl	45
46	30% (v/v) PEG-200	CAPS pH 10.5	0.2 M (NH ₄) ₂ SO ₄	46
47	50% (v/v) PEG-200	HEPES pH 7.5	none	47
48	50% (v/v) PEG-200	phosphate-citrate pH 4.2	0.2 M NaCl	48

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