

**Wizard I random sparse matrix crystallization screen - technical sheet**

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

	<u>crystallant</u>	<u>buffer (0.1 M)</u>	<u>salt (0.2 M)</u>	
1	20% (w/v) PEG-8000	CHES pH 9.5	none	1
2	10% (v/v) 2-propanol	HEPES pH 7.5	NaCl	2
3	15% (v/v) ethanol	CHES pH 9.5	none	3
4	35% (v/v) 2-methyl-2,4-pentanediol	imidazole pH 8.0	MgCl <sub>2</sub>	4
5	30% (v/v) PEG-400	CAPS pH 10.5	none	5
6	20% (w/v) PEG-3000	citrate pH 5.5	none	6
7	10% (w/v) PEG-8000	MES pH 6.0	Zn(OAc) <sub>2</sub>	7
8	2.0 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	citrate pH 5.5	none	8
9	1.0 M (NH <sub>4</sub> ) <sub>2</sub> HPO <sub>4</sub>	acetate pH 4.5	none	9
10	20% (w/v) PEG-2000 MME	Tris pH 7.0	none	10
11	20% (v/v) 1,4-butanediol	MES pH 6.0	Li <sub>2</sub> SO <sub>4</sub>	11
12	20% (w/v) PEG-1000	imidazole pH 8.0	Ca(OAc) <sub>2</sub>	12
13	1.26 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	cacodylate pH 6.5	none	13
14	1.0 M sodium citrate	cacodylate pH 6.5	none	14
15	10% (w/v) PEG-3000	imidazole pH 8.0	Li <sub>2</sub> SO <sub>4</sub>	15
16	2.5 M NaCl	Na/K phosphate pH 6.2	none	16
17	30% (w/v) PEG-8000	acetate pH 4.5	Li <sub>2</sub> SO <sub>4</sub>	17
18	1.0 M K/Na tartrate	imidazole pH 8.0	NaCl	18
19	20% (w/v) PEG-1000	Tris pH 7.0	none	19
20	0.4 M NaH <sub>2</sub> PO <sub>4</sub> /1.6 M K <sub>2</sub> HPO <sub>4</sub>	imidazole pH 8.0	NaCl	20
21	20% (w/v) PEG-8000	HEPES pH 7.5	none	21
22	10% (v/v) 2-propanol	Tris pH 8.5	none	22
23	15% (v/v) ethanol	imidazole pH 8.0	MgCl <sub>2</sub>	23
24	35% (v/v) 2-methyl-2,4-pentanediol	Tris pH 7.0	NaCl	24
25	30% (v/v) PEG-400	Tris pH 8.5	MgCl <sub>2</sub>	25
26	10% (w/v) PEG-3000	CHES pH 9.5	none	26
27	1.2 M NaH <sub>2</sub> PO <sub>4</sub> /0.8 M K <sub>2</sub> HPO <sub>4</sub>	CAPS pH 10.5	Li <sub>2</sub> SO <sub>4</sub>	27
28	20% (w/v) PEG-3000	HEPES pH 7.5	NaCl	28
29	10% (w/v) PEG-8000	CHES pH 9.5	NaCl	29
30	1.26 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	acetate pH 4.5	NaCl	30
31	20% (w/v) PEG-8000	phosphate-citrate pH 4.2	NaCl	31
32	10% (w/v) PEG-3000	Na/K phosphate pH 6.2	none	32
33	2.0 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	CAPS pH 10.5	Li <sub>2</sub> SO <sub>4</sub>	33
34	1.0 M (NH <sub>4</sub> ) <sub>2</sub> HPO <sub>4</sub>	imidazole pH 8.0	none	34
35	20% (v/v) 1,4-butanediol	acetate pH 4.5	none	35
36	1.0 M sodium citrate	imidazole pH 8.0	none	36
37	2.5 M NaCl	imidazole pH 8.0	none	37
38	1.0 M K/Na tartrate	CHES pH 9.5	Li <sub>2</sub> SO <sub>4</sub>	38
39	20% (w/v) PEG-1000	phosphate-citrate pH 4.2	Li <sub>2</sub> SO <sub>4</sub>	39
40	10% (v/v) 2-propanol	MES pH 6.0	Ca(OAc) <sub>2</sub>	40
41	30% (w/v) PEG-3000	CHES pH 9.5	none	41
42	15% (v/v) ethanol	Tris pH 7.0	none	42
43	35% (v/v) 2-methyl-2,4-pentanediol	Na/K phosphate pH 6.2	none	43
44	30% (v/v) PEG-400	acetate pH 4.5	Ca(OAc) <sub>2</sub>	44
45	20% (w/v) PEG-3000	acetate pH 4.5	none	45
46	10% (w/v) PEG-8000	imidazole pH 8.0	Ca(OAc) <sub>2</sub>	46
47	1.26 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	Tris pH 8.5	Li <sub>2</sub> SO <sub>4</sub>	47
48	20% (w/v) PEG-1000	acetate pH 4.5	Zn(OAc) <sub>2</sub>	48

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

## Wizard II random sparse matrix crystallization screen - technical sheet

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

	<u>crystallant</u>	<u>buffer (0.1 M)</u>	<u>salt (0.2 M)</u>	
1	10% (w/v) PEG-3000	acetate pH 4.5	Zn(OAc) <sub>2</sub>	1
2	35% (v/v) 2-methyl-2,4-pentanediol	MES pH 6.0	Li <sub>2</sub> SO <sub>4</sub>	2
3	20% (w/v) PEG-8000	Tris pH 8.5	MgCl <sub>2</sub>	3
4	2.0 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	cacodylate pH 6.5	NaCl	4
5	20% (v/v) 1,4-butanediol	HEPES pH 7.5	NaCl	5
6	10% (v/v) 2-propanol	phosphate-citrate pH 4.2	Li <sub>2</sub> SO <sub>4</sub>	6
7	30% (w/v) PEG-3000	Tris pH 7.0	NaCl	7
8	10% (w/v) PEG-8000	Na/K phosphate pH 6.2	NaCl	8
9	2.0 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	phosphate-citrate pH 4.2	none	9
10	1.0 M (NH <sub>4</sub> ) <sub>2</sub> HPO <sub>4</sub>	Tris pH 8.5	none	10
11	10% (v/v) 2-propanol	cacodylate pH 6.5	Zn(OAc) <sub>2</sub>	11
12	30% (v/v) PEG-400	cacodylate pH 6.5	Li <sub>2</sub> SO <sub>4</sub>	12
13	15% (v/v) ethanol	citrate pH 5.5	Li <sub>2</sub> SO <sub>4</sub>	13
14	20% (w/v) PEG-1000	Na/K phosphate pH 6.2	NaCl	14
15	1.26 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	HEPES pH 7.5	none	15
16	1.0 M sodium citrate	CHES pH 9.5	none	16
17	2.5 M NaCl	Tris pH 7.0	MgCl <sub>2</sub>	17
18	20% (w/v) PEG-3000	Tris pH 7.0	Ca(OAc) <sub>2</sub>	18
19	1.6 M NaH <sub>2</sub> PO <sub>4</sub> /0.4 M K <sub>2</sub> HPO <sub>4</sub>	phosphate-citrate pH 4.2	none	19
20	15% (v/v) ethanol	MES pH 6.0	Zn(OAc) <sub>2</sub>	20
21	35% (v/v) 2-methyl-2,4-pentanediol	acetate pH 4.5	none	21
22	10% (v/v) 2-propanol	imidazole pH 8.0	none	22
23	15% (v/v) ethanol	HEPES pH 7.5	MgCl <sub>2</sub>	23
24	30% (w/v) PEG-8000	imidazole pH 8.0	NaCl	24
25	35% (v/v) 2-methyl-2,4-pentanediol	HEPES pH 7.5	NaCl	25
26	30% (v/v) PEG-400	CHES pH 9.5	none	26
27	10% (w/v) PEG-3000	cacodylate pH 6.5	MgCl <sub>2</sub>	27
28	20% (w/v) PEG-8000	MES pH 6.0	Ca(OAc) <sub>2</sub>	28
29	1.26 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	CHES pH 9.5	NaCl	29
30	20% (v/v) 1,4-butanediol	imidazole pH 8.0	Zn(OAc) <sub>2</sub>	30
31	1.0 M sodium citrate	Tris pH 7.0	NaCl	31
32	20% (w/v) PEG-1000	Tris pH 8.5	none	32
33	1.0 M (NH <sub>4</sub> ) <sub>2</sub> HPO <sub>4</sub>	citrate pH 5.5	NaCl	33
34	10% (w/v) PEG-8000	imidazole pH 8.0	none	34
35	0.8 M NaH <sub>2</sub> PO <sub>4</sub> /1.2 M K <sub>2</sub> HPO <sub>4</sub>	acetate pH 4.5	none	35
36	10% (w/v) PEG-3000	phosphate-citrate pH 4.2	NaCl	36
37	1.0 M K/Na tartrate	Tris pH 7.0	Li <sub>2</sub> SO <sub>4</sub>	37
38	2.5 M NaCl	acetate pH 4.5	Li <sub>2</sub> SO <sub>4</sub>	38
39	20% (w/v) PEG-8000	CAPS pH 10.5	NaCl	39
40	20% (w/v) PEG-3000	imidazole pH 8.0	Zn(OAc) <sub>2</sub>	40
41	2.0 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	Tris pH 7.0	Li <sub>2</sub> SO <sub>4</sub>	41
42	30% (v/v) PEG-400	HEPES pH 7.5	NaCl	42
43	10% (w/v) PEG-8000	Tris pH 7.0	MgCl <sub>2</sub>	43
44	20% (w/v) PEG-1000	cacodylate pH 6.5	MgCl <sub>2</sub>	44
45	1.26 M (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>	MES pH 6.0	none	45
46	1.0 M (NH <sub>4</sub> ) <sub>2</sub> HPO <sub>4</sub>	imidazole pH 8.0	NaCl	46
47	2.5 M NaCl	imidazole pH 8.0	Zn(OAc) <sub>2</sub>	47
48	1.0 M K/Na tartrate	MES pH 6.0	none	48

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