

**PEG-Anomalous™ PEG 1K 48-Salt crystallization screen - technical sheet**

**Formulations: (Patent Pending)**

	<u>PEG-1K</u>	<u>anomalous salt</u>
1	30%(w/v) PEG 1000	0.2M ammonium bromide
2	30%(w/v) PEG 1000	0.2M ammonium iodide
3	30%(w/v) PEG 1000	0.2M ammonium selenate
4	30%(w/v) PEG 1000	0.2M barium acetate
5	30%(w/v) PEG 1000	0.2M barium bromide, dihydrate
6	30%(w/v) PEG 1000	0.2M barium chloride, dihydrate
7	30%(w/v) PEG 1000	0.2M barium iodide
8	30%(w/v) PEG 1000	0.2M cadmium acetate
9	30%(w/v) PEG 1000	0.2M cadmium bromide, tetrahydrate
10	30%(w/v) PEG 1000	0.2M cadmium iodide
11	30%(w/v) PEG 1000	0.2M cadmium nitrate
12	30%(w/v) PEG 1000	0.2M cadmium sulfate, hydrate
13	30%(w/v) PEG 1000	0.2M cesium acetate
14	30%(w/v) PEG 1000	0.2M cesium bromide
15	30%(w/v) PEG 1000	0.2M cesium chloride
16	30%(w/v) PEG 1000	0.2M cesium fluoride
17	30%(w/v) PEG 1000	0.2M cesium formate
18	30%(w/v) PEG 1000	0.2M cesium iodide
19	30%(w/v) PEG 1000	0.2M cesium nitrate
20	30%(w/v) PEG 1000	0.2M cesium sulfate
21	30%(w/v) PEG 1000	0.2M gadolinium acetate, hydrate
22	30%(w/v) PEG 1000	0.2M gadolinium bromide, hydrate
23	30%(w/v) PEG 1000	0.2M gadolinium chloride, hydrate
24	30%(w/v) PEG 1000	0.2M gadolinium nitrate
25	30%(w/v) PEG 1000	0.2M holmium bromide
26	30%(w/v) PEG 1000	0.2M holmium chloride
27	30%(w/v) PEG 1000	0.2M holmium nitrate
28	30%(w/v) PEG 1000	0.2M lanthanum acetate, hydrate
29	30%(w/v) PEG 1000	0.2M Lanthanum bromide
30	30%(w/v) PEG 1000	0.2M lanthanum chloride
31	30%(w/v) PEG 1000	0.2M Lanthanum nitrate
32	30%(w/v) PEG 1000	0.2M lithium bromide
33	30%(w/v) PEG 1000	0.2M lithium iodide
34	30%(w/v) PEG 1000	0.2M magnesium bromide
35	30%(w/v) PEG 1000	0.2M potassium bromide
36	30%(w/v) PEG 1000	0.2M potassium iodide
37	30%(w/v) PEG 1000	0.2M potassium selenate
38	30%(w/v) PEG 1000	0.2M samarium bromide, hexahydrate
39	30%(w/v) PEG 1000	0.2M samarium chloride, hexahydrate
40	30%(w/v) PEG 1000	0.2M samarium nitrate
41	30%(w/v) PEG 1000	0.2M sodium bromide
42	30%(w/v) PEG 1000	0.2M sodium iodide
43	30%(w/v) PEG 1000	0.2M sodium selenate
44	30%(w/v) PEG 1000	0.2M sodium tungstate
45	30%(w/v) PEG 1000	0.2M Strontium nitrate
46	30%(w/v) PEG 1000	0.2M ytterbium chloride
47	30%(w/v) PEG 1000	0.2M Yttrium nitrate
48	30%(w/v) PEG 1000	0.2M zinc bromide

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

**PEG-Anomalous™ PEG 4K 48-Salt crystallization screen - technical sheet****Formulations: (Patent Pending)**

	<u>PEG-4K</u>	<u>anomalous salt</u>
1	20%(w/v) PEG 4000	0.2M ammonium bromide
2	20%(w/v) PEG 4000	0.2M ammonium iodide
3	20%(w/v) PEG 4000	0.2M ammonium selenate
4	20%(w/v) PEG 4000	0.2M barium acetate
5	20%(w/v) PEG 4000	0.2M barium bromide, dihydrate
6	20%(w/v) PEG 4000	0.2M barium chloride, dihydrate
7	20%(w/v) PEG 4000	0.2M barium iodide
8	20%(w/v) PEG 4000	0.2M cadmium acetate
9	20%(w/v) PEG 4000	0.2M cadmium bromide, tetrahydrate
10	20%(w/v) PEG 4000	0.2M cadmium iodide
11	20%(w/v) PEG 4000	0.2M cadmium nitrate
12	20%(w/v) PEG 4000	0.2M cadmium sulfate, hydrate
13	20%(w/v) PEG 4000	0.2M cesium acetate
14	20%(w/v) PEG 4000	0.2M cesium bromide
15	20%(w/v) PEG 4000	0.2M cesium chloride
16	20%(w/v) PEG 4000	0.2M cesium fluoride
17	20%(w/v) PEG 4000	0.2M cesium formate
18	20%(w/v) PEG 4000	0.2M cesium iodide
19	20%(w/v) PEG 4000	0.2M cesium nitrate
20	20%(w/v) PEG 4000	0.2M cesium sulfate
21	20%(w/v) PEG 4000	0.2M gadolinium acetate, hydrate
22	20%(w/v) PEG 4000	0.2M gadolinium bromide, hydrate
23	20%(w/v) PEG 4000	0.2M gadolinium chloride, hydrate
24	20%(w/v) PEG 4000	0.2M gadolinium nitrate
25	20%(w/v) PEG 4000	0.2M holmium bromide
26	20%(w/v) PEG 4000	0.2M holmium chloride
27	20%(w/v) PEG 4000	0.2M holmium nitrate
28	20%(w/v) PEG 4000	0.2M lanthanum acetate, hydrate
29	20%(w/v) PEG 4000	0.2M Lanthanum bromide
30	20%(w/v) PEG 4000	0.2M lanthanum chloride
31	20%(w/v) PEG 4000	0.2M Lanthanum nitrate
32	20%(w/v) PEG 4000	0.2M lithium bromide
33	20%(w/v) PEG 4000	0.2M lithium iodide
34	20%(w/v) PEG 4000	0.2M magnesium bromide
35	20%(w/v) PEG 4000	0.2M potassium bromide
36	20%(w/v) PEG 4000	0.2M potassium iodide
37	20%(w/v) PEG 4000	0.2M potassium selenate
38	20%(w/v) PEG 4000	0.2M samarium bromide, hexahydrate
39	20%(w/v) PEG 4000	0.2M samarium chloride, hexahydrate
40	20%(w/v) PEG 4000	0.2M samarium nitrate
41	20%(w/v) PEG 4000	0.2M sodium bromide
42	20%(w/v) PEG 4000	0.2M sodium iodide
43	20%(w/v) PEG 4000	0.2M sodium selenate
44	20%(w/v) PEG 4000	0.2M sodium tungstate
45	20%(w/v) PEG 4000	0.2M Strontium nitrate
46	20%(w/v) PEG 4000	0.2M ytterbium chloride
47	20%(w/v) PEG 4000	0.2M Yttrium nitrate
48	20%(w/v) PEG 4000	0.2M zinc bromide

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](http://www.emeraldbiosystems.com)

**PEG-Anomalous™ PEG 8K 48-Salt crystallization screen - technical sheet**

**Formulations: (Patent Pending)**

	<u>PEG-8K</u>	<u>anomalous salt</u>
1	20%(w/v) PEG 8000	0.2M ammonium bromide
2	20%(w/v) PEG 8000	0.2M ammonium iodide
3	20%(w/v) PEG 8000	0.2M ammonium selenate
4	20%(w/v) PEG 8000	0.2M barium acetate
5	20%(w/v) PEG 8000	0.2M barium bromide, dihydrate
6	20%(w/v) PEG 8000	0.2M barium chloride, dihydrate
7	20%(w/v) PEG 8000	0.2M barium iodide
8	20%(w/v) PEG 8000	0.2M cadmium acetate
9	20%(w/v) PEG 8000	0.2M cadmium bromide, tetrahydrate
10	20%(w/v) PEG 8000	0.2M cadmium iodide
11	20%(w/v) PEG 8000	0.2M cadmium nitrate
12	20%(w/v) PEG 8000	0.2M cadmium sulfate, hydrate
13	20%(w/v) PEG 8000	0.2M cesium acetate
14	20%(w/v) PEG 8000	0.2M cesium bromide
15	20%(w/v) PEG 8000	0.2M cesium chloride
16	20%(w/v) PEG 8000	0.2M cesium fluoride
17	20%(w/v) PEG 8000	0.2M cesium formate
18	20%(w/v) PEG 8000	0.2M cesium iodide
19	20%(w/v) PEG 8000	0.2M cesium nitrate
20	20%(w/v) PEG 8000	0.2M cesium sulfate
21	20%(w/v) PEG 8000	0.2M gadolinium acetate, hydrate
22	20%(w/v) PEG 8000	0.2M gadolinium bromide, hydrate
23	20%(w/v) PEG 8000	0.2M gadolinium chloride, hydrate
24	20%(w/v) PEG 8000	0.2M gadolinium nitrate
25	20%(w/v) PEG 8000	0.2M holmium bromide
26	20%(w/v) PEG 8000	0.2M holmium chloride
27	20%(w/v) PEG 8000	0.2M holmium nitrate
28	20%(w/v) PEG 8000	0.2M lanthanum acetate, hydrate
29	20%(w/v) PEG 8000	0.2M Lanthanum bromide
30	20%(w/v) PEG 8000	0.2M lanthanum chloride
31	20%(w/v) PEG 8000	0.2M Lanthanum nitrate
32	20%(w/v) PEG 8000	0.2M lithium bromide
33	20%(w/v) PEG 8000	0.2M lithium iodide
34	20%(w/v) PEG 8000	0.2M magnesium bromide
35	20%(w/v) PEG 8000	0.2M potassium bromide
36	20%(w/v) PEG 8000	0.2M potassium iodide
37	20%(w/v) PEG 8000	0.2M potassium selenate
38	20%(w/v) PEG 8000	0.2M samarium bromide, hexahydrate
39	20%(w/v) PEG 8000	0.2M samarium chloride, hexahydrate
40	20%(w/v) PEG 8000	0.2M samarium nitrate
41	20%(w/v) PEG 8000	0.2M sodium bromide
42	20%(w/v) PEG 8000	0.2M sodium iodide
43	20%(w/v) PEG 8000	0.2M sodium selenate
44	20%(w/v) PEG 8000	0.2M sodium tungstate
45	20%(w/v) PEG 8000	0.2M Strontium nitrate
46	20%(w/v) PEG 8000	0.2M ytterbium chloride
47	20%(w/v) PEG 8000	0.2M Yttrium nitrate
48	20%(w/v) PEG 8000	0.2M zinc bromide

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

**PEG-Anomalous™ PEG 10K 48-Salt crystallization screen - technical sheet**

**Formulations: (Patent Pending)**

	<u>PEG-10K</u>	<u>anomalous salt</u>
1	10%(w/v) PEG 10,000	0.2M ammonium bromide
2	10%(w/v) PEG 10,000	0.2M ammonium iodide
3	10%(w/v) PEG 10,000	0.2M ammonium selenate
4	10%(w/v) PEG 10,000	0.2M barium acetate
5	10%(w/v) PEG 10,000	0.2M barium bromide, dihydrate
6	10%(w/v) PEG 10,000	0.2M barium chloride, dihydrate
7	10%(w/v) PEG 10,000	0.2M barium iodide
8	10%(w/v) PEG 10,000	0.2M cadmium acetate
9	10%(w/v) PEG 10,000	0.2M cadmium bromide, tetrahydrate
10	10%(w/v) PEG 10,000	0.2M cadmium iodide
11	10%(w/v) PEG 10,000	0.2M cadmium nitrate
12	10%(w/v) PEG 10,000	0.2M cadmium sulfate, hydrate
13	10%(w/v) PEG 10,000	0.2M cesium acetate
14	10%(w/v) PEG 10,000	0.2M cesium bromide
15	10%(w/v) PEG 10,000	0.2M cesium chloride
16	10%(w/v) PEG 10,000	0.2M cesium fluoride
17	10%(w/v) PEG 10,000	0.2M cesium formate
18	10%(w/v) PEG 10,000	0.2M cesium iodide
19	10%(w/v) PEG 10,000	0.2M cesium nitrate
20	10%(w/v) PEG 10,000	0.2M cesium sulfate
21	10%(w/v) PEG 10,000	0.2M gadolinium acetate, hydrate
22	10%(w/v) PEG 10,000	0.2M gadolinium bromide, hydrate
23	10%(w/v) PEG 10,000	0.2M gadolinium chloride, hydrate
24	10%(w/v) PEG 10,000	0.2M gadolinium nitrate
25	10%(w/v) PEG 10,000	0.2M holmium bromide
26	10%(w/v) PEG 10,000	0.2M holmium chloride
27	10%(w/v) PEG 10,000	0.2M holmium nitrate
28	10%(w/v) PEG 10,000	0.2M lanthanum acetate, hydrate
29	10%(w/v) PEG 10,000	0.2M Lanthanum bromide
30	10%(w/v) PEG 10,000	0.2M lanthanum chloride
31	10%(w/v) PEG 10,000	0.2M Lanthanum nitrate
32	10%(w/v) PEG 10,000	0.2M lithium bromide
33	10%(w/v) PEG 10,000	0.2M lithium iodide
34	10%(w/v) PEG 10,000	0.2M magnesium bromide
35	10%(w/v) PEG 10,000	0.2M potassium bromide
36	10%(w/v) PEG 10,000	0.2M potassium iodide
37	10%(w/v) PEG 10,000	0.2M potassium selenate
38	10%(w/v) PEG 10,000	0.2M samarium bromide, hexahydrate
39	10%(w/v) PEG 10,000	0.2M samarium chloride, hexahydrate
40	10%(w/v) PEG 10,000	0.2M samarium nitrate
41	10%(w/v) PEG 10,000	0.2M sodium bromide
42	10%(w/v) PEG 10,000	0.2M sodium iodide
43	10%(w/v) PEG 10,000	0.2M sodium selenate
44	10%(w/v) PEG 10,000	0.2M sodium tungstate
45	10%(w/v) PEG 10,000	0.2M Strontium nitrate
46	10%(w/v) PEG 10,000	0.2M ytterbium chloride
47	10%(w/v) PEG 10,000	0.2M Yttrium nitrate
48	10%(w/v) PEG 10,000	0.2M zinc bromide

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