

**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