

SmartScreen – PurePEGS



The PurePEGS screen features mixtures of purified PEGs, from 300 MW to 8,000 MW, in each condition at a total concentration of 22.5 % (1) combined with salts and buffers giving the most crystallization hits spanning a wide pH range (1.1 - 9.8).

Features:

- **Ultra pure PEGs – USP grade**
- **Optimized for diffusive mixing; capillary crystallization**
- **Polydisperse PEGs in each condition**
- **pH measured and recorded for final solution**
- **Better control of excipients in the PEGs**

Papers for additional information:

1. J. Newman *et al.*, *Acta Crystallogr D Biol Crystallogr* **61**, 1426 (2005).
2. R. Page *et al.*, *Acta Crystallogr D Biol Crystallogr* **59**, 1028 (2003).
3. J. W. Wooh, R. D. Kidd, J. L. Martin, B. Kobe, *Acta Crystallogr D Biol Crystallogr* **59**, 769 (2003).
4. J. T. Bukrinsky, J. C. N. Poulsen, *Journal of Applied Crystallography* **34**, 533 (2001).

MLY#	Position	Composition	Final pH
1	A1	0.3 M Ammonium chloride , 0.1 M Citric acid , 22.5% PurePEGS PEG 0.3-8kD	4
2	A2	0.3 M Ammonium formate , 0.1 M Potassium nitrate , 22.5% PurePEGS PEG 0.3-8kD	6.6
3	B1	0.3 M Ammonium nitrate , 0.1 M Sodium citrate tribasic dihydrate , 22.5% PurePEGS PEG 0.3-8kD	5.9
4	B2	0.3 M Ammonium phosphate monobasic, 0.1 M MES monohydrate , 22.5% PurePEGS PEG 0.3-8kD	5.7
5	C1	0.3 M Ammonium sulfate , 0.1 M BIS-TRIS , 22.5% PurePEGS PEG 0.3-8kD	6.8
6	C2	0.3 M Di-ammonium tartrate , 0.1 M Sodium acetate trihydrate , 22.5% PurePEGS PEG 0.3-8kD	5.8
7	D1	0.3 M Cadmium chloride , 0.1 M Sodium acetate trihydrate , 22.5% PurePEGS PEG 0.3-8kD	3.7
8	D2	0.3 M Cadmium sulfate , 0.1 M Imidazole , 22.5% PurePEGS PEG 0.3-8kD	4.6
9	E1	0.3 M Caesium chloride, 0.1 M HEPES , 22.5% PurePEGS PEG 0.3-8kD	7.6
10	E2	0.3 M Calcium acetate, 0.1 M BICINE , 22.5% PurePEGS PEG 0.3-8kD	6.4
11	F1	0.3 M Calcium chloride, 0.1 M BIS-TRIS propane , 22.5% PurePEGS PEG 0.3-8kD	8.5
12	F2	0.3 M Cobalt(II) chloride hexahydrate, 0.1 M Ammonium citrate, 22.5% PurePEGS PEG 0.3-8kD	3.2
13	G1	0.3 M Di-ammonium hydrogen citrate , 0.1 M Magnesium formate, 22.5% PurePEGS PEG	4.9

		0.3-8kD	
14	G2	0.3 M DL- malic acid, 0.1 M Sodium cacodylate trihydrate, 22.5% PurePEGs PEG 0.3-8kD	2.9
15	H1	0.3 M Iron(III) chloride hexahydrate, 0.1 M Potassium nitrate, 22.5% PurePEGs PEG 0.3-8kD	1.1
16	H2	0.3 M Lithium chloride, 0.1 M Potassium formate, 22.5% PurePEGs PEG 0.3-8kD	6.3
17	A6	0.3 M Lithium sulfate monohydrate, 0.1 M CHES, 22.5% PurePEGs PEG 0.3-8kD	9.5
18	A7	0.3 M Magnesium acetate tetrahydrate, 0.1 M Citric acid , 22.5% PurePEGs PEG 0.3-8kD	4.7
19	B6	0.3 M Magnesium chloride hexahydrate, 0.1 M Sodium cacodylate trihydrate, 22.5% PurePEGs PEG 0.3-8kD	5.8
20	B7	0.3 M Magnesium nitrate hexahydrate, 0.1 M HEPES , 22.5% PurePEGs PEG 0.3-8kD	7.4
21	C6	0.3 M Magnesium sulfate heptahydrate, 0.1 M Potassium formate, 22.5% PurePEGs PEG 0.3-8kD	6.1
22	C7	0.3 M Nickel(II) chloride hexahydrate, 0.1 M MES monohydrate , 22.5% PurePEGs PEG 0.3-8kD	5.7
23	D6	0.3 M Potassium bromide, 0.1 M BIS-TRIS , 22.5% PurePEGs PEG 0.3-8kD	6.6
24	D7	0.3 M Potassium phosphate monobasic, 0.1 M Tris , 22.5% PurePEGs PEG 0.3-8kD	6.1
25	E6	0.3 M Potassium sodium tartrate tetrahydrate, 0.1 M BICINE , 22.5% PurePEGs PEG 0.3-8kD	8.8
26	E7	0.3 M Potassium tartrate, 0.1 M Imidazole , 22.5% PurePEGs PEG 0.3-8kD	7.3
27	F6	0.3 M Potassium thiocyanate, 0.1 M Ammonium citrate, 22.5% PurePEGs PEG 0.3-8kD	5.8
28	F7	0.3 M Sodium bromide, 0.1 M Sodium cacodylate trihydrate, 22.5% PurePEGs PEG 0.3-8kD	6.5
29	G6	0.3 M Sodium chloride, 0.1 M Tris , 22.5% PurePEGs PEG 0.3-8kD	8.1
30	G7	0.3 M Sodium fluoride, 0.1 M CHES, 22.5% PurePEGs PEG 0.3-8kD	9.6
31	H6	0.3 M Sodium formate, 0.1 M Citric acid , 22.5% PurePEGs PEG 0.3-8kD	4.5
32	H7	0.3 M Sodium iodide, 0.1 M Sodium acetate trihydrate , 22.5% PurePEGs PEG 0.3-8kD	4.7
33	A11	0.3 M Sodium malonate, 0.1 M Sodium citrate tribasic dihydrate , 22.5% PurePEGs PEG 0.3-8kD	6.8
34	A12	0.3 M Sodium nitrate, 0.1 M MES monohydrate , 22.5% PurePEGs PEG 0.3-8kD	6.1
35	B11	0.3 M Sodium phosphate monobasic monohydrate, 0.1 M BIS-TRIS , 22.5% PurePEGs PEG 0.3-8kD	5.8
36	B12	0.3 M Sodium sulfate, 0.1 M BIS-TRIS propane , 22.5% PurePEGs PEG 0.3-8kD	9.3
37	C11	0.3 M Sodium thiocyanate , 0.1 M HEPES , 22.5% PurePEGs PEG 0.3-8kD	7.8
38	C12	0.3 M Succinic acid , 0.1 M Imidazole , 22.5% PurePEGs PEG 0.3-8kD	3.4
39	D11	0.3 M Tri-methylamine N-oxide, 0.1 M BICINE , 22.5% PurePEGs PEG 0.3-8kD	8.6
40	D12	0.3 M Tri-potassium citrate , 0.1 M BIS-TRIS propane , 22.5% PurePEGs PEG 0.3-8kD	9.8
41	E11	10 % Jeffamine M-600, 0.1 M HEPES , 22.5% PurePEGs PEG 0.3-8kD	7.5
42	E12	0.3 M Zinc chloride, 0.1 M Magnesium formate, 22.5% PurePEGs PEG 0.3-8kD	5.7
43	F11	0.3 M Zinc sulfate heptahydrate, 0.1 M Sodium citrate tribasic dihydrate , 22.5% PurePEGs PEG 0.3-8kD	4
44	F12	0.1 M CHES, 22.5% PurePEGs PEG 0.3-8kD	9.9
45	G11	0.1 M Imidazole , 22.5% PurePEGs PEG 0.3-8kD	6.6
46	G12	0.1 M Magnesium formate, 22.5% PurePEGs PEG 0.3-8kD	5.9
47	H11	0.1 M Sodium acetate trihydrate , 22.5% PurePEGs PEG 0.3-8kD	4.7
48	H12	0.1 M Tris , 22.5% PurePEGs PEG 0.3-8kD	8.1