



Cat.# EBS-BWZ-EXT

Wizard™IV random sparse matrix crystallization screen - technical sheet

Formulations:

	crystallant	buffer (0.1 M)	additive(s)	additive(s)	
1 (A1)	20%(v/v) Glycerol		40mM Potassium phosphate	16%(w/v) PEG 8000	1 (A1)
2 (A2)	15%(v/v) Ethanol	Tris	8.0 100mM Sodium chloride	5%(v/v) MPD	2 (A2)
3 (A3)	40%(v/v) Ethanol	Phosphate-citrate	4.2	5%(w/v) PEG 1000	3 (A3)
4 (A4)	200mM Ammonium sulfate	BisTris	5.5		4 (A4)
5 (A5)	2M Ammonium sulfate	Acetate	5.5	2%(v/v) PEG 400	5 (A5)
6 (A6)	800mM Ammonium Sulfate	Citrate	4.0		6 (A6)
7 (A7)	2M Lithium sulfate	Acetate	4.5	100mM Magnesium sulfate	7 (A7)
8 (A8)	2M Lithium sulfate	Tris	8.5	2%(v/v) PEG 400	8 (A8)
9 (A9)	2M Lithium sulfate	Acetate	5.5	100mM Magnesium sulfate	9 (A9)
10 (A10)	50%(v/v) PEG 200	Cacodylate	6.5	200mM Magnesium chloride	10 (A10)
11 (A11)	40%(v/v) PEG 300	Cacodylate	6.5	200mM Calcium acetate	11 (A11)
12 (A12)	30%(v/v) Jeffamine M600 pH 7.0	HEPES	7.0		12 (A12)
13 (B1)	800mM Succinic Acid pH 7.0				13 (B1)
14 (B2)	40%(v/v) PEG 400	Tris	8.5	200mM Lithium sulfate	14 (B2)
15 (B3)	50%(v/v) PEG 400	Acetate	4.5	200mM Lithium sulfate	15 (B3)
16 (B4)	15%(v/v) PEG 550MME	MES	6.5		16 (B4)
17 (B5)	25%(w/v) PEG 1500	SPG Buffer/NaOH	5.5		17 (B5)
18 (B6)	25%(w/v) PEG 1500	SPG Buffer/NaOH	8.5		18 (B6)
19 (B7)	25%(w/v) PEG 1500	MMT Buffer/NaOH	6.5		19 (B7)
20 (B8)	25%(w/v) PEG 1500	MMT Buffer/NaOH	9.0		20 (B8)
21 (B9)	25%(w/v) PEG 1500	MIB Buffer/HCl	5.0		21 (B9)
22 (B10)	25%(w/v) PEG 1500	PCB Buffer/NaOH	7.0		22 (B10)
23 (B11)	12%(w/v) PEG 1500	Acetate	5.5	2500mM Sodium chloride	23 (B11)
24 (B12)	2400mM Sodium Malonate			1.5%(v/v) MPD	24 (B12)
25 (C1)	30%(w/v) PEG 2000MME			150mM Potassium bromide	25 (C1)
26 (C2)	10%(w/v) PEG 2000MME	Sodium acetate	5.5	200mM Ammonium sulfate	26 (C2)
27 (C3)	20%(w/v) PEG 2000MME	Tris	8.5	200mM Trimethylamine n-oxide	27 (C3)
28 (C4)	20%(w/v) PEG 3350	BisTris Propane	6.5	200mM Sodium fluoride	28 (C4)
29 (C5)	20%(w/v) PEG 3350	Citrate	4.0	200mM Sodium citrate	29 (C5)
30 (C6)	20%(w/v) PEG 3350	BisTris Propane	8.5	200mM Sodium malonate	30 (C6)
31 (C7)	20%(w/v) Polyacrylic acid 5100	HEPES	7.0	20mM Magnesium chloride	31 (C7)
32 (C8)	2100mM DL Malic acid pH 7.0				32 (C8)
33 (C9)	800mM Potassium phosphate (dibasic)	HEPES	7.5	800mM Sodium Phosphate	33 (C9)
34 (C10)	20%(w/v) PEG 6000	MES	6.0	200mM Ammonium chloride	34 (C10)
35 (C11)	20%(w/v) PEG 6000	HEPES	7.0	200mM Sodium chloride	35 (C11)
36 (C12)	20%(w/v) PEG 6000	Tris	8.0	200mM Lithium chloride	36 (C12)
37 (D1)	20%(w/v) Polyvinylpyrrolidone K15	Tris	8.5	100mM Cobalt chloride	37 (D1)
38 (D2)	50%(v/v) Ethylene glycol	Tris	8.5	200mM Magnesium chloride	38 (D2)
39 (D3)	20%(w/v) PEG 8000	Imidazole	6.5	3%(v/v) MPD	39 (D3)
40 (D4)	20%(w/v) PEG 8000	Tris	8.5	100mM Magnesium chloride	40 (D4)
41 (D5)	20%(w/v) PEG 8000	HEPES	7.5	200mM Ammonium sulfate	41 (D5)
42 (D6)	30%(v/v) MPD	Acetate	4.5	25%(w/v) PEG 1500	42 (D6)
43 (D7)	30%(v/v) MPD	Imidazole	6.5	10%(w/v) PEG 3350	43 (D7)
44 (D8)	30%(v/v) MPD	Tris	8.5	500mM Sodium chloride	44 (D8)
45 (D9)	40%(v/v) Isopropanol	Imidazole	6.5	15%(w/v) PEG 8000	45 (D9)
46 (D10)	30%(v/v) Isopropanol	Tris	8.5	30%(w/v) PEG 3350	46 (D10)
47 (D11)	17%(w/v) PEG 10000	BisTris	5.5	100mM Ammonium Acetate	47 (D11)
48 (D12)	15%(w/v) PEG 20000	HEPES	7.0		48 (D12)

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

© 2009 Emerald BioSystems, Inc., (888) 780-8535, FAX (206) 780-8549, www.emeraldbiosystems.com