



**Wizard™ III random sparse matrix crystallization screen - technical sheet**

**Formulations:**

|    | <u>crystallant</u>                | <u>buffer</u>       | <u>pH</u> | <u>salt/additive #1</u>              | <u>additive #2</u> |
|----|-----------------------------------|---------------------|-----------|--------------------------------------|--------------------|
| 1  | 20%(w/v) PEG 3350                 |                     |           | .2M Ammonium citrate (dibasic)       |                    |
| 2  | 30%(v/v) MPD                      | .1M Sodium acetate  | 4.6       | .02M Calcium chloride                |                    |
| 3  | 20%(w/v) PEG 3350                 |                     |           | .2M Magnesium formate                |                    |
| 4  | 20%(w/v) PEG 3350                 |                     |           | .2M Ammonium formate                 |                    |
| 5  | 20%(w/v) PEG 3350                 |                     |           | .2M Ammonium chloride                |                    |
| 6  | 20%(w/v) PEG 3350                 |                     |           | .2M Potassium formate                |                    |
| 7  | 50%(w/v) MPD                      | .1M Tris            | 8.5       | .2M Ammonium phosphate (monobasic)   |                    |
| 8  | 20%(w/v) PEG 3350                 |                     |           | .2M Potassium nitrate                |                    |
| 9  | 0.8M Ammonium sulfate             | .1M Citric acid     | 4.0       |                                      |                    |
| 10 | 20%(w/v) PEG 3350                 |                     |           | .2M Sodium thiocyanate               |                    |
| 11 | 20%(w/v) PEG 6000                 | .1M Bicine          | 9.0       |                                      |                    |
| 12 | 10%(w/v) PEG 8000                 | .1M HEPES           | 7.5       | 8%(v/v) Ethylene glycol              |                    |
| 13 | 8%(w/v) PEG 4000                  | .1M Sodium acetate  | 4.6       |                                      |                    |
| 14 | 20%(w/v) PEG 6000                 | .1M Citric acid     | 5.0       |                                      |                    |
| 15 | 1.6M Sodium citrate               |                     |           |                                      |                    |
| 16 | 20%(w/v) PEG 3350                 |                     |           | .2M Potassium citrate tribasic       |                    |
| 17 | 20%(w/v) PEG 4000                 | .1M Citrate         | 5.5       | 10%(v/v) 2-propanol                  |                    |
| 18 | 20%(w/v) PEG 6000                 | .1M Citric acid     | 4.0       | 1M Lithium chloride                  |                    |
| 19 | 20%(w/v) PEG 3350                 |                     |           | .2M Ammonium nitrate                 |                    |
| 20 | 10%(w/v) PEG 6000                 | .1M HEPES           | 7.0       |                                      |                    |
| 21 | .8M Potassium phosphate (dibasic) | .1M HEPES           | 7.5       | .8M Sodium phosphate (monobasic)     |                    |
| 22 | 20%(v/v) Ethanol                  | .1M Tris            | 8.5       |                                      |                    |
| 23 | 10%(w/v) PEG 20,000               | .1M Bicine          | 9.0       | 2%(v/v) Dioxane                      |                    |
| 24 | 2M Ammonium sulfate               | .1M Sodium acetate  | 4.6       |                                      |                    |
| 25 | 10%(w/v) PEG 1000                 |                     |           | 10%(w/v) PEG 8000                    |                    |
| 26 | 24%(w/v) PEG 1500                 |                     |           | 20%(v/v) Glycerol                    |                    |
| 27 | 30%(v/v) PEG 400                  | .1M HEPES           | 7.5       | .2M Magnesium chloride               |                    |
| 28 | 70%(v/v) MPD                      | .1M HEPES           | 7.5       |                                      |                    |
| 29 | 40%(v/v) MPD                      | .1M Tris            | 8.0       |                                      |                    |
| 30 | 25.5%(w/v) PEG 4000               |                     |           | .17M Ammonium sulfate                | 15%(v/v) Glycerol  |
| 31 | 14%(v/v) 2-propanol               | .07M Sodium acetate | 4.6       | .14M Calcium chloride                | 30%(v/v) Glycerol  |
| 32 | 16%(w/v) PEG 8000                 |                     |           | .04M Potassium phosphate (monobasic) | 20%(v/v) Glycerol  |
| 33 | 1.6M Magnesium sulfate            | .1M MES             | 6.5       |                                      |                    |
| 34 | 10%(w/v) PEG 6000                 | .1M Bicine          | 9.0       |                                      |                    |
| 35 | 14.4%(w/v) PEG 8000               | .08M Cacodylate     | 6.5       | .16M Calcium acetate                 | 20%(v/v) Glycerol  |
| 36 | 30%(v/v) Jeffamine M-600          | .1M MES             | 6.5       | .05M Cesium chloride                 |                    |
| 37 | 3.2M Ammonium sulfate             | .1M Citric acid     | 5.0       |                                      |                    |
| 38 | 15%(w/v) PEG 10,000               | .1M Citrate         | 5.5       | 2%(v/v) Dioxane                      |                    |
| 39 | 20%(v/v) Jeffamine M-600          | .1M HEPES           | 7.5       |                                      |                    |
| 40 | 10%(v/v) MPD                      | .1M Bicine          | 9.0       |                                      |                    |
| 41 | 28%(v/v) PEG 400                  | .1M HEPES           | 7.5       | .2M Calcium chloride                 |                    |
| 42 | 30%(w/v) PEG 4000                 | .1M Tris            | 8.5       | .2M Lithium sulfate                  |                    |
| 43 | 30%(w/v) PEG 8000                 |                     |           | .2M Ammonium sulfate                 |                    |
| 44 | 30%(w/v) PEG 5000 MME             | .1M Tris            | 8.0       | .2M Lithium sulfate                  |                    |
| 45 | 1.5M Ammonium sulfate             | .1M Tris            | 8.5       |                                      | 12%(v/v) Glycerol  |
| 46 | 50%(v/v) MPD                      | .1M Tris            | 8.5       | .2M Ammonium chloride                |                    |
| 47 | 30%(w/v) PEG 5000 MME             | .1M MES             | 6.5       | .2M Ammonium sulfate                 |                    |
| 48 | 20%(w/v) PEG 10,000               | .1M HEPES           | 7.5       |                                      |                    |

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



Cat.# EBS-WIZ-EXT

**Wizard™IV random sparse matrix crystallization screen - technical sheet**

**Formulations:**

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

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