

**PEG-Anomalous™ PEG 10K 48-Salt crystallization screen - technical sheet****Formulations: (Patent Pending)**

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