

## Ozma PEG 1K 48-Salt crystallization screen - technical sheet

### Formulations:

	<u>Precipitant</u>		<u>Salt</u>	
<b>1</b>	30%(w/v) PEG 1000	200 mM	Ammonium acetate	<b>1</b>
<b>2</b>	30%(w/v) PEG 1000	200 mM	Ammonium chloride	<b>2</b>
<b>3</b>	30%(w/v) PEG 1000	200 mM	Ammonium fluoride	<b>3</b>
<b>4</b>	30%(w/v) PEG 1000	200 mM	Ammonium formate	<b>4</b>
<b>5</b>	30%(w/v) PEG 1000	200 mM	Ammonium iodide	<b>5</b>
<b>6</b>	30%(w/v) PEG 1000	200 mM	Ammonium nitrate	<b>6</b>
<b>7</b>	30%(w/v) PEG 1000	200 mM	Ammonium phosphate (monobasic)	<b>7</b>
<b>8</b>	30%(w/v) PEG 1000	200 mM	Ammonium phosphate (dibasic)	<b>8</b>
<b>9</b>	30%(w/v) PEG 1000	200 mM	Ammonium sulfate	<b>9</b>
<b>10</b>	30%(w/v) PEG 1000	200 mM	Ammonium sulfite	<b>10</b>
<b>11</b>	30%(w/v) PEG 1000	200 mM	Calcium acetate	<b>11</b>
<b>12</b>	30%(w/v) PEG 1000	200 mM	Calcium chloride	<b>12</b>
<b>13</b>	30%(w/v) PEG 1000	200 mM	Diammonium tartrate	<b>13</b>
<b>14</b>	30%(w/v) PEG 1000	200 mM	Formic acid potassium salt	<b>14</b>
<b>15</b>	30%(w/v) PEG 1000	200 mM	Lithium acetate	<b>15</b>
<b>16</b>	30%(w/v) PEG 1000	200 mM	Lithium chloride	<b>16</b>
<b>17</b>	30%(w/v) PEG 1000	200 mM	Lithium citrate	<b>17</b>
<b>18</b>	30%(w/v) PEG 1000	200 mM	Lithium nitrate	<b>18</b>
<b>19</b>	30%(w/v) PEG 1000	200 mM	Lithium sulfate	<b>19</b>
<b>20</b>	30%(w/v) PEG 1000	200 mM	Magnesium acetate	<b>20</b>
<b>21</b>	30%(w/v) PEG 1000	200 mM	Magnesium chloride	<b>21</b>
<b>22</b>	30%(w/v) PEG 1000	200 mM	Magnesium formate	<b>22</b>
<b>23</b>	30%(w/v) PEG 1000	200 mM	Magnesium nitrate	<b>23</b>
<b>24</b>	30%(w/v) PEG 1000	200 mM	Magnesium sulfate	<b>24</b>
<b>25</b>	30%(w/v) PEG 1000	200 mM	Potassium acetate	<b>25</b>
<b>26</b>	30%(w/v) PEG 1000	200 mM	Potassium chloride	<b>26</b>
<b>27</b>	30%(w/v) PEG 1000	200 mM	Potassium fluoride	<b>27</b>
<b>28</b>	30%(w/v) PEG 1000	200 mM	Potassium iodide	<b>28</b>
<b>29</b>	30%(w/v) PEG 1000	200 mM	Potassium nitrate	<b>29</b>
<b>30</b>	30%(w/v) PEG 1000	50 mM	Potassium phosphate (dibasic)	<b>30</b>
<b>31</b>	30%(w/v) PEG 1000	200 mM	Potassium phosphate (monobasic)	<b>31</b>
<b>32</b>	30%(w/v) PEG 1000	200 mM	Potassium sulfate	<b>32</b>
<b>33</b>	30%(w/v) PEG 1000	200 mM	Potassium thiocyanate	<b>33</b>
<b>34</b>	30%(w/v) PEG 1000	200 mM	Potassium/Sodium tartrate	<b>34</b>
<b>35</b>	30%(w/v) PEG 1000	200 mM	Sodium acetate	<b>35</b>
<b>36</b>	30%(w/v) PEG 1000	150 mM	Sodium chloride	<b>36</b>
<b>37</b>	30%(w/v) PEG 1000	200 mM	Sodium citrate	<b>37</b>
<b>38</b>	30%(w/v) PEG 1000	200 mM	Sodium fluoride	<b>38</b>
<b>39</b>	30%(w/v) PEG 1000	200 mM	Sodium formate	<b>39</b>
<b>40</b>	30%(w/v) PEG 1000	200 mM	Sodium iodide	<b>40</b>
<b>41</b>	30%(w/v) PEG 1000	200 mM	Sodium isothiocyanate	<b>41</b>
<b>42</b>	30%(w/v) PEG 1000	200 mM	Sodium nitrate	<b>42</b>
<b>43</b>	30%(w/v) PEG 1000	200 mM	Sodium phosphate (dibasic)	<b>43</b>
<b>44</b>	30%(w/v) PEG 1000	200 mM	Sodium phosphate (monobasic)	<b>44</b>
<b>45</b>	30%(w/v) PEG 1000	200 mM	Sodium sulfate	<b>45</b>
<b>46</b>	30%(w/v) PEG 1000	200 mM	Sodium tartrate (dibasic)	<b>46</b>
<b>47</b>	30%(w/v) PEG 1000	200 mM	Tripotassium citrate	<b>47</b>
<b>48</b>	30%(w/v) PEG 1000	200 mM	Zinc acetate	<b>48</b>

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)

## Ozma PEG 4K 48-Salt crystallization screen - technical sheet

### Formulations:

	<u>Precipitant</u>		<u>Salt</u>	
<b>1</b>	20%(w/v) PEG 4000	200 mM	Ammonium acetate	<b>1</b>
<b>2</b>	20%(w/v) PEG 4000	200 mM	Ammonium chloride	<b>2</b>
<b>3</b>	20%(w/v) PEG 4000	200 mM	Ammonium fluoride	<b>3</b>
<b>4</b>	20%(w/v) PEG 4000	200 mM	Ammonium formate	<b>4</b>
<b>5</b>	20%(w/v) PEG 4000	200 mM	Ammonium iodide	<b>5</b>
<b>6</b>	20%(w/v) PEG 4000	200 mM	Ammonium nitrate	<b>6</b>
<b>7</b>	20%(w/v) PEG 4000	200 mM	Ammonium phosphate (monobasic)	<b>7</b>
<b>8</b>	20%(w/v) PEG 4000	200 mM	Ammonium phosphate (dibasic)	<b>8</b>
<b>9</b>	20%(w/v) PEG 4000	200 mM	Ammonium sulfate	<b>9</b>
<b>10</b>	20%(w/v) PEG 4000	200 mM	Ammonium sulfite	<b>10</b>
<b>11</b>	20%(w/v) PEG 4000	200 mM	Calcium acetate	<b>11</b>
<b>12</b>	20%(w/v) PEG 4000	50 mM	Calcium chloride	<b>12</b>
<b>13</b>	20%(w/v) PEG 4000	200 mM	Diammonium tartrate	<b>13</b>
<b>14</b>	20%(w/v) PEG 4000	200 mM	Formic acid potassium salt	<b>14</b>
<b>15</b>	20%(w/v) PEG 4000	200 mM	Lithium acetate	<b>15</b>
<b>16</b>	20%(w/v) PEG 4000	200 mM	Lithium chloride	<b>16</b>
<b>17</b>	20%(w/v) PEG 4000	200 mM	Lithium citrate	<b>17</b>
<b>18</b>	20%(w/v) PEG 4000	200 mM	Lithium nitrate	<b>18</b>
<b>19</b>	20%(w/v) PEG 4000	200 mM	Lithium sulfate	<b>19</b>
<b>20</b>	20%(w/v) PEG 4000	200 mM	Magnesium acetate	<b>20</b>
<b>21</b>	20%(w/v) PEG 4000	50 mM	Magnesium chloride	<b>21</b>
<b>22</b>	20%(w/v) PEG 4000	200 mM	Magnesium formate	<b>22</b>
<b>23</b>	20%(w/v) PEG 4000	200 mM	Magnesium nitrate	<b>23</b>
<b>24</b>	20%(w/v) PEG 4000	200 mM	Magnesium sulfate	<b>24</b>
<b>25</b>	20%(w/v) PEG 4000	200 mM	Potassium acetate	<b>25</b>
<b>26</b>	20%(w/v) PEG 4000	100 mM	Potassium chloride	<b>26</b>
<b>27</b>	20%(w/v) PEG 4000	200 mM	Potassium fluoride	<b>27</b>
<b>28</b>	20%(w/v) PEG 4000	200 mM	Potassium iodide	<b>28</b>
<b>29</b>	20%(w/v) PEG 4000	200 mM	Potassium nitrate	<b>29</b>
<b>30</b>	20%(w/v) PEG 4000	200 mM	Potassium phosphate (dibasic)	<b>30</b>
<b>31</b>	20%(w/v) PEG 4000	100 mM	Potassium phosphate (monobasic)	<b>31</b>
<b>32</b>	20%(w/v) PEG 4000	100 mM	Potassium sulfate	<b>32</b>
<b>33</b>	20%(w/v) PEG 4000	200 mM	Potassium thiocyanate	<b>33</b>
<b>34</b>	20%(w/v) PEG 4000	200 mM	Potassium/Sodium tartrate	<b>34</b>
<b>35</b>	20%(w/v) PEG 4000	100 mM	Sodium acetate	<b>35</b>
<b>36</b>	20%(w/v) PEG 4000	200 mM	Sodium chloride	<b>36</b>
<b>37</b>	20%(w/v) PEG 4000	50 mM	Sodium citrate	<b>37</b>
<b>38</b>	20%(w/v) PEG 4000	200 mM	Sodium fluoride	<b>38</b>
<b>39</b>	20%(w/v) PEG 4000	200 mM	Sodium formate	<b>39</b>
<b>40</b>	20%(w/v) PEG 4000	200 mM	Sodium iodide	<b>40</b>
<b>41</b>	20%(w/v) PEG 4000	200 mM	Sodium isothiocyanate	<b>41</b>
<b>42</b>	20%(w/v) PEG 4000	200 mM	Sodium nitrate	<b>42</b>
<b>43</b>	20%(w/v) PEG 4000	200 mM	Sodium phosphate (dibasic)	<b>43</b>
<b>44</b>	20%(w/v) PEG 4000	50 mM	Sodium phosphate (monobasic)	<b>44</b>
<b>45</b>	20%(w/v) PEG 4000	200 mM	Sodium sulfate	<b>45</b>
<b>46</b>	20%(w/v) PEG 4000	200 mM	Sodium tartrate (dibasic)	<b>46</b>
<b>47</b>	20%(w/v) PEG 4000	100 mM	Tripotassium citrate	<b>47</b>
<b>48</b>	20%(w/v) PEG 4000	200 mM	Zinc acetate	<b>48</b>

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)

## **Ozma PEG 8K 48-Salt crystallization screen - technical sheet**

### **Formulations:**

	<b><u>precipitant</u></b>		<b><u>Salt</u></b>	
<b>1</b>	20%(w/v) PEG 8000	200 mM	Ammonium acetate	<b>1</b>
<b>2</b>	20%(w/v) PEG 8000	200 mM	Ammonium chloride	<b>2</b>
<b>3</b>	20%(w/v) PEG 8000	200 mM	Ammonium fluoride	<b>3</b>
<b>4</b>	20%(w/v) PEG 8000	200 mM	Ammonium formate	<b>4</b>
<b>5</b>	20%(w/v) PEG 8000	200 mM	Ammonium iodide	<b>5</b>
<b>6</b>	20%(w/v) PEG 8000	200 mM	Ammonium nitrate	<b>6</b>
<b>7</b>	20%(w/v) PEG 8000	200 mM	Ammonium phosphate (monobasic)	<b>7</b>
<b>8</b>	20%(w/v) PEG 8000	200 mM	Ammonium phosphate (dibasic)	<b>8</b>
<b>9</b>	20%(w/v) PEG 8000	200 mM	Ammonium sulfate	<b>9</b>
<b>10</b>	20%(w/v) PEG 8000	200 mM	Ammonium sulfite	<b>10</b>
<b>11</b>	20%(w/v) PEG 8000	200 mM	Calcium acetate	<b>11</b>
<b>12</b>	20%(w/v) PEG 8000	50 mM	Calcium chloride	<b>12</b>
<b>13</b>	20%(w/v) PEG 8000	200 mM	Diammonium tartrate	<b>13</b>
<b>14</b>	20%(w/v) PEG 8000	200 mM	Formic acid potassium salt	<b>14</b>
<b>15</b>	20%(w/v) PEG 8000	200 mM	Lithium acetate	<b>15</b>
<b>16</b>	20%(w/v) PEG 8000	200 mM	Lithium chloride	<b>16</b>
<b>17</b>	20%(w/v) PEG 8000	200 mM	Lithium citrate	<b>17</b>
<b>18</b>	20%(w/v) PEG 8000	200 mM	Lithium nitrate	<b>18</b>
<b>19</b>	20%(w/v) PEG 8000	200 mM	Lithium sulfate	<b>19</b>
<b>20</b>	20%(w/v) PEG 8000	200 mM	Magnesium acetate	<b>20</b>
<b>21</b>	20%(w/v) PEG 8000	200 mM	Magnesium chloride	<b>21</b>
<b>22</b>	20%(w/v) PEG 8000	200 mM	Magnesium formate	<b>22</b>
<b>23</b>	20%(w/v) PEG 8000	200 mM	Magnesium nitrate	<b>23</b>
<b>24</b>	20%(w/v) PEG 8000	200 mM	Magnesium sulfate	<b>24</b>
<b>25</b>	20%(w/v) PEG 8000	200 mM	Potassium acetate	<b>25</b>
<b>26</b>	20%(w/v) PEG 8000	200 mM	Potassium chloride	<b>26</b>
<b>27</b>	20%(w/v) PEG 8000	200 mM	Potassium fluoride	<b>27</b>
<b>28</b>	20%(w/v) PEG 8000	200 mM	Potassium iodide	<b>28</b>
<b>29</b>	20%(w/v) PEG 8000	200 mM	Potassium nitrate	<b>29</b>
<b>30</b>	20%(w/v) PEG 8000	200 mM	Potassium phosphate (dibasic)	<b>30</b>
<b>31</b>	20%(w/v) PEG 8000	100 mM	Potassium phosphate (monobasic)	<b>31</b>
<b>32</b>	20%(w/v) PEG 8000	200 mM	Potassium sulfate	<b>32</b>
<b>33</b>	20%(w/v) PEG 8000	200 mM	Potassium thiocyanate	<b>33</b>
<b>34</b>	20%(w/v) PEG 8000	200 mM	Potassium/Sodium tartrate	<b>34</b>
<b>35</b>	20%(w/v) PEG 8000	50 mM	Sodium acetate	<b>35</b>
<b>36</b>	20%(w/v) PEG 8000	200 mM	Sodium chloride	<b>36</b>
<b>37</b>	20%(w/v) PEG 8000	100 mM	Sodium citrate	<b>37</b>
<b>38</b>	20%(w/v) PEG 8000	200 mM	Sodium fluoride	<b>38</b>
<b>39</b>	20%(w/v) PEG 8000	200 mM	Sodium formate	<b>39</b>
<b>40</b>	20%(w/v) PEG 8000	200 mM	Sodium iodide	<b>40</b>
<b>41</b>	20%(w/v) PEG 8000	200 mM	Sodium isothiocyanate	<b>41</b>
<b>42</b>	20%(w/v) PEG 8000	200 mM	Sodium nitrate	<b>42</b>
<b>43</b>	20%(w/v) PEG 8000	200 mM	Sodium phosphate (dibasic)	<b>43</b>
<b>44</b>	20%(w/v) PEG 8000	100 mM	Sodium phosphate (monobasic)	<b>44</b>
<b>45</b>	20%(w/v) PEG 8000	200 mM	Sodium sulfate	<b>45</b>
<b>46</b>	20%(w/v) PEG 8000	200 mM	Sodium tartrate (dibasic)	<b>46</b>
<b>47</b>	20%(w/v) PEG 8000	200 mM	Tripotassium citrate	<b>47</b>
<b>48</b>	20%(w/v) PEG 8000	200 mM	Zinc acetate	<b>48</b>

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)

## Ozma PEG 10K 48-Salt crystallization screen - technical sheet

### Formulations:

	<u>precipitant</u>		<u>salt</u>	
<b>1</b>	10%(w/v) PEG 10,000	200 mM	Ammonium acetate	<b>1</b>
<b>2</b>	10%(w/v) PEG 10,000	200 mM	Ammonium chloride	<b>2</b>
<b>3</b>	10%(w/v) PEG 10,000	200 mM	Ammonium fluoride	<b>3</b>
<b>4</b>	10%(w/v) PEG 10,000	200 mM	Ammonium formate	<b>4</b>
<b>5</b>	10%(w/v) PEG 10,000	200 mM	Ammonium iodide	<b>5</b>
<b>6</b>	10%(w/v) PEG 10,000	200 mM	Ammonium nitrate	<b>6</b>
<b>7</b>	10%(w/v) PEG 10,000	200 mM	Ammonium phosphate (monobasic)	<b>7</b>
<b>8</b>	10%(w/v) PEG 10,000	200 mM	Ammonium phosphate (dibasic)	<b>8</b>
<b>9</b>	10%(w/v) PEG 10,000	200 mM	Ammonium sulfate	<b>9</b>
<b>10</b>	10%(w/v) PEG 10,000	200 mM	Ammonium sulfite	<b>10</b>
<b>11</b>	10%(w/v) PEG 10,000	200 mM	Calcium acetate	<b>11</b>
<b>12</b>	10%(w/v) PEG 10,000	200 mM	Calcium chloride	<b>12</b>
<b>13</b>	10%(w/v) PEG 10,000	200 mM	Diammonium tartrate	<b>13</b>
<b>14</b>	10%(w/v) PEG 10,000	200 mM	Formic acid potassium salt	<b>14</b>
<b>15</b>	10%(w/v) PEG 10,000	200 mM	Lithium acetate	<b>15</b>
<b>16</b>	10%(w/v) PEG 10,000	200 mM	Lithium chloride	<b>16</b>
<b>17</b>	10%(w/v) PEG 10,000	200 mM	Lithium citrate	<b>17</b>
<b>18</b>	10%(w/v) PEG 10,000	200 mM	Lithium nitrate	<b>18</b>
<b>19</b>	10%(w/v) PEG 10,000	200 mM	Lithium sulfate	<b>19</b>
<b>20</b>	10%(w/v) PEG 10,000	200 mM	Magnesium acetate	<b>20</b>
<b>21</b>	10%(w/v) PEG 10,000	200 mM	Magnesium chloride	<b>21</b>
<b>22</b>	10%(w/v) PEG 10,000	200 mM	Magnesium formate	<b>22</b>
<b>23</b>	10%(w/v) PEG 10,000	200 mM	Magnesium nitrate	<b>23</b>
<b>24</b>	10%(w/v) PEG 10,000	200 mM	Magnesium sulfate	<b>24</b>
<b>25</b>	10%(w/v) PEG 10,000	200 mM	Potassium acetate	<b>25</b>
<b>26</b>	10%(w/v) PEG 10,000	200 mM	Potassium chloride	<b>26</b>
<b>27</b>	10%(w/v) PEG 10,000	200 mM	Potassium fluoride	<b>27</b>
<b>28</b>	10%(w/v) PEG 10,000	200 mM	Potassium iodide	<b>28</b>
<b>29</b>	10%(w/v) PEG 10,000	200 mM	Potassium nitrate	<b>29</b>
<b>30</b>	10%(w/v) PEG 10,000	200 mM	Potassium phosphate (dibasic)	<b>30</b>
<b>31</b>	10%(w/v) PEG 10,000	50 mM	Potassium phosphate (monobasic)	<b>31</b>
<b>32</b>	10%(w/v) PEG 10,000	200 mM	Potassium sulfate	<b>32</b>
<b>33</b>	10%(w/v) PEG 10,000	200 mM	Potassium thiocyanate	<b>33</b>
<b>34</b>	10%(w/v) PEG 10,000	200 mM	Potassium/Sodium tartrate	<b>34</b>
<b>35</b>	10%(w/v) PEG 10,000	200 mM	Sodium acetate	<b>35</b>
<b>36</b>	10%(w/v) PEG 10,000	150 mM	Sodium chloride	<b>36</b>
<b>37</b>	10%(w/v) PEG 10,000	200 mM	Sodium citrate	<b>37</b>
<b>38</b>	10%(w/v) PEG 10,000	200 mM	Sodium fluoride	<b>38</b>
<b>39</b>	10%(w/v) PEG 10,000	200 mM	Sodium formate	<b>39</b>
<b>40</b>	10%(w/v) PEG 10,000	200 mM	Sodium iodide	<b>40</b>
<b>41</b>	10%(w/v) PEG 10,000	200 mM	Sodium isothiocyanate	<b>41</b>
<b>42</b>	10%(w/v) PEG 10,000	200 mM	Sodium nitrate	<b>42</b>
<b>43</b>	10%(w/v) PEG 10,000	200 mM	Sodium phosphate (dibasic)	<b>43</b>
<b>44</b>	10%(w/v) PEG 10,000	200 mM	Sodium phosphate (monobasic)	<b>44</b>
<b>45</b>	10%(w/v) PEG 10,000	200 mM	Sodium sulfate	<b>45</b>
<b>46</b>	10%(w/v) PEG 10,000	200 mM	Sodium tartrate (dibasic)	<b>46</b>
<b>47</b>	10%(w/v) PEG 10,000	200 mM	Tripotassium citrate	<b>47</b>
<b>48</b>	10%(w/v) PEG 10,000	200 mM	Zinc acetate	<b>48</b>

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)