



**A Division of Prolume Ltd.**

www.Nanolight.com info@prolume.com Tel: 928-367-1200

## **Instructions: #3692 Water-soluble *Prolume Purple* for *in vitro* use**

### **Content:**

231.8 µg water soluble Prolume Purple (MW: 463.53) per vial  
resulting in a 50 µM Prolume Purple solution  
if dissolved in 10 ml of an aqueous buffer

### **Storage:**

Please store vacuum sealed vials at -80°C, for shorter periods of time at -20°C. Keep dark.  
Dissolved Prolume Purple should be used immediately.

### **Usage:**

1. Warm tube to room temperature.
2. For a 50 µM solution add 10 ml of an aqueous buffer (e.g. PBS, TBS).  
For a 100 µM solution add only 5 ml.
3. Vortex until completely dissolved. Use immediately in a luminometer assay.

Please note that the stability of the compound is dependent on the pH of the added buffer. Buffers with a pH greater than 6.5 will cause a faster degradation (oxidation) of the Coelenterazine compound.