

Instructions for CAT#3032 Water-soluble Coelenterazine for *in vitro* use

Content:

212 µg native Coelenterazine/vial with water-solubilizing additives resulting in a 50 µM Coelenterazine-Solution if dissolved in 10 ml of an aqueous buffer

Storage:

Please store unopened vials at -80°C, for shorter periods of time at -20°C. Keep dark and dry. Dissolved Coelenterazine should be used immediately.

Usage:

1. Warm vial to room temperature.
2. For a 50 µM solution add 10 ml of an aqueous buffer (e.g. PBS).
3. Wait 30 sec. and invert vial a few times until completely dissolved. Use immediately in a luminometer assay.

Application sample:

