



A Division of Prolume Ltd.

www.nanolight.com

info@prolume.com

Tel: 928-367-1200

Instructions: #3702 Water-soluble v-Coelenterazine for *in vitro* use

Content:

224 µg v-Coelenterazine/tube with water solubilizing additives
resulting in a 50 µM v-Coelenterazine-Solution
if dissolved in 10 ml of an aqueous buffer

Storage:

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

Usage:

1. Warm tube to room temperature. (Do NOT use a heat block.)
2. For a 50 µM solution add 10 ml of an aqueous buffer (e.g. PBS), for a more concentrated solution add less.
3. Wait for 30 sec. and invert a few times until completely dissolved. Use immediately in a luminometer assay.

Note: v-Coelenterazine is in general more instable than native Coelenterazine. Once dissolved in an aqueous solution use promptly. Buffers with a pH lower than 7.0 will increase the long-term stability.