

## Manual for h-Coelenterazine

Cat.# 301 h-Coelenterazine CAS# 50909-86-9 MW 407.46

**Alternative names:** Renilla Luciferin, 2-Deoxycoelenterazine, Coelenterazine-h

**General Notes:** h-Coelenterazine (h-CTZ) is the dehydroxy derivative of native Coelenterazine and can be used to increase the initial luminous intensity of Aequorin up to 20-fold, thus becoming a more sensitive  $\text{Ca}^{2+}$  sensor.

Dr O. Shimomura (Biochem. J. 1989 Aug 1;261(3); 913-920) was able to show that h-Coelenterazine produced 10 times more initial light intensity but the total number of photons produced was 82% overall. Essentially the initial light output was compressed and 10 times the flash intensity was noted over a moderately shorter rise time.

h-Coelenterazine is also used as substrate for Renilla Luciferase, creating a lower initial intensity but a longer lasting kinetic.

h-Coelenterazine should not be used with Gaussia Luciferase, it will require native Coelenterazine (Cat.#303).

**Storage and Storage Life:** It is best stored as completely dry powder under Argon in air-tight O-ring plastic tubes at  $-20^{\circ}\text{C}$  or for longer storage at  $-80^{\circ}\text{C}$ , protected from light. Oxygen and moisture will lead to auto-oxidation of CTZ over time, reducing its overall activity.

**Dissolving h-Coelenterazine:** We recommend using our specifically developed NanoFuel Solvent (Cat. #399) for maximum solubility and shelf-life. Adding  $500\ \mu\text{l}$  to  $500\ \mu\text{g}$  of lyophilized h-Coelenterazine will result in a  $1\ \text{mg/ml}$  solution that can be stored at  $-20^{\circ}\text{C}$  or below for at least one year without any notable degradation.

As an alternative you may use Methanol to dissolve h-Coelenterazine. To prevent oxidation, it is recommended to acidify and degas the alcohol prior to addition.

**Dilution and luminescent assaying:** We recommend to prepare the working solution fresh every time before a luminometer assay. In general, a  $100\ \mu\text{M}$  h-Coelenterazine solution will work with most assays.

Use  $407.5\ \mu\text{l}$  of the  $1\ \text{mg/ml}$  stock solution and dilute in  $10\ \text{ml}$  of your buffer of choice (e.g. PBS) to get a  $100\ \mu\text{M}$  solution. All solutions should be at room temperature. When performing a plate reader assay and you want to compare relative light units (RLU) results between the first well and last well in a 96 plate, please note that h-CTZ will continuously oxidize over time in aqueous solutions (see graph, performed with two widely used buffers).

