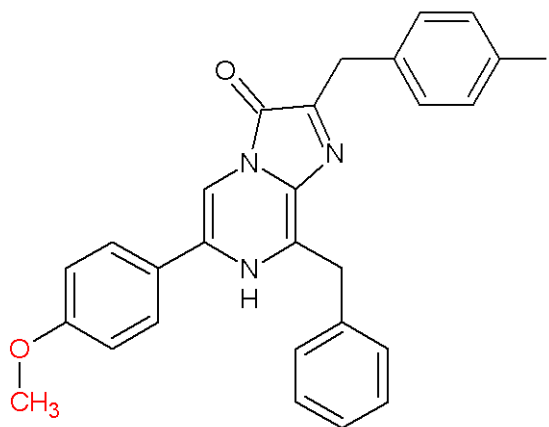


## HOW TO USE “Prolume Purple IV”

Cat. #363    Methoxy-COELENTERAZINE-Iodine    patent pending    MW 547.39



**General Notes:** Methoxy Coelenterazine-I (Me-O-CTZ-I) called “Prolume Purple IV” is a new synthetic analogue of Coelenterazine-I with an additional **methoxy group**. This compound was developed by Nanolight™ Technology and has not tested yet with all available luciferases.

**Storage and Shelf Life:** It is best stored as completely DRY powder under argon in air-tight O-ring plastic tubes at -20°C or for longer storage at -70°C, protected from light.

**Usage:** It is always best to make FRESH SOLUTIONS immediately before luminometer assays or experiments.

1. Dissolve lyophilized **Prolume Purple IV** in **NanoFuel Solvent** as a 1 mg/ml solution (Ethanol won't dissolve Me-O-CTZ-I).
2. Use this stock solution to make an aqueous solution in PBS or TBS (e.g. 50 µM for luminometer assays equal to 273.7 µl (1mg/ml) in 10 ml PBS).
3. Store dissolved **Prolume Purple IV** at -80°C, do not store the aqueous working solution (it will oxidize over time). **Prolume Purple IV** has the same stability in aqueous solutions like any other Coelenterazine analogue.