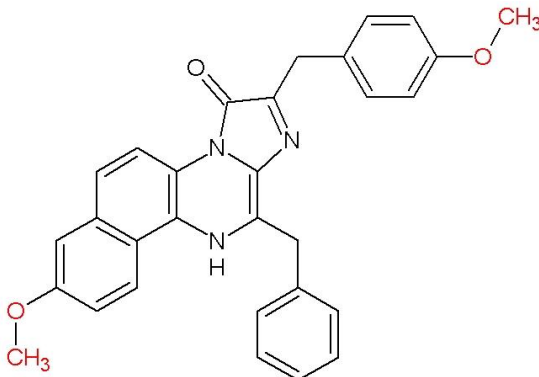


HOW TO USE “Prolume Purple V”

Cat. #361 Methoxy-v-COELENTERAZINE-Methoxy patent pending MW 475.19



General Notes: Methoxy-v-Coelenterazine-Methoxy (Me-O-v-CTZ-O-Me) called “Prolume Purple V” is a new synthetic analogue of v-Coelenterazine with **two** 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 V** in **NanoFuel Solvent** as a 1 mg/ml solution (Ethanol won't dissolve Me-O-v-CTZ-O-Me).
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 V** at -80°C, do not store the aqueous working solution (it will oxidize over time). **Prolume Purple V** has the same stability in aqueous solutions like any other Coelenterazine analogue.