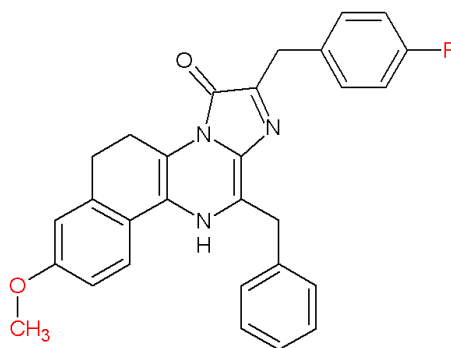


HOW TO USE “Prolume Purple III”

Cat. #365 Methoxy e-COELENTERAZINE-F patent pending MW 465.52



General Notes: Methoxy e-Coelenterazine-F (Me-O-e-CTZ-F) called “**Prolume Purple III**” is a new synthetic analogue of e-Coelenterazine-F with an additional **methoxy group**. This compound was developed by Nanolight™ Technology to work with **Green Renilla Luciferase (Green RLuc)** to emit light at **~410 nm**.

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.

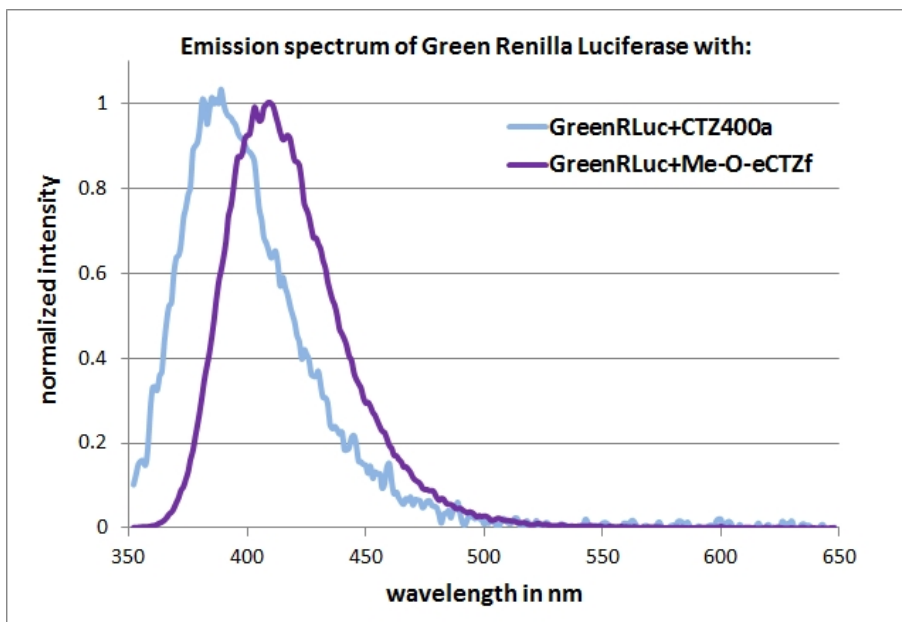
Usage: It is always best to make FRESH SOLUTIONS immediately before luminometer assays or experiments.

1. Dissolve lyophilized **Prolume Purple III** in **NanoFuel Solvent** as a 1 mg/ml solution (Ethanol won't dissolve Me-O-eCTZF).
2. Use this stock solution to make an aqueous solution in PBS or TBS (e.g. 50 μM for luminometer assays equal to 232.8 μl (1mg/ml) in 10 ml PBS).
3. Store dissolved **Prolume Purple III** at -80°C, do not store the aqueous working solution (it will oxidize over time). **Prolume Purple III** has the same stability in aqueous solutions like any other Coelenterazine analogue.

HOW TO USE “Prolume Purple III”

Properties of Nanolight™ Prolume Purple III Cat. #365

A. Comparison of emission spectra of e-CTZ and “Prolume Purple” with Green *Renilla muelleri* Luciferase



B. Activity of CTZ analogues with Green *Renilla muelleri* Luciferase

