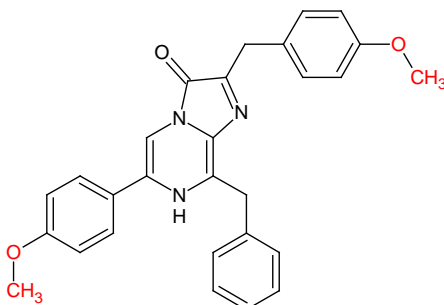


## Certificate of Analysis

### CAT# 367 “Prolume Purple II” Methoxy-Coelenterazine-Methoxy

Date of Test: 09 October 2019  
 Compound ID: Methoxy-Coelenterazine-Methoxy  
 CAS#: Not Listed, patent pending  
 IUPAC Name: Me-O-CTZ-O-Me or “Prolume Purple II”  
 Batch Number: J2913 (11606-045)  
 Manufacturer: Prolume Ltd  
 Molecular Formula: C<sub>28</sub>H<sub>25</sub>N<sub>3</sub>O<sub>3</sub>  
 MW: 451.52 g/mol

Structure:



Test Items	Specification	Results	Test method
<b>Appearance</b>	Pale Yellow Solid	Conforms	Visual
<b>Purity</b>	>95%	98.6%	HPLC
<b>Identification</b>	NMR is consistent with structure	Conforms	NMR
	Mass Spec is consistent with structure	Conforms	MS
<b>Storage</b>	Store dry and dark at -20°C or for long term storage at -80°C.		