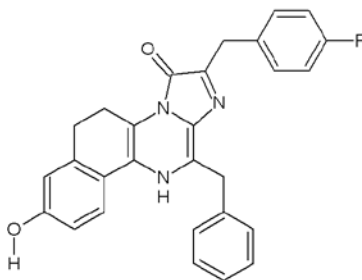


## Certificate of Analysis

### CAT# 346 e-Coelenterazine-F\*

Date of Test: 13 December 2019  
 Compound ID: eCTZ-F  
 CAS#: Not registered  
 IUPAC Name: e-Coelenterazine-Fluoride  
 Batch Number: 996-03611-150  
 Manufacturer: Prolume Ltd  
 Molecular Formula: C<sub>28</sub>H<sub>22</sub>FN<sub>3</sub>O<sub>2</sub>  
 MW: 451.49 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Orange Solid	Conforms	Visual
Purity	>95%	95.6%	HPLC
Identification	Spectrum is consistent with structure	Conforms	By NMR
	Emission at 413nm and 475nm	Conforms	By Activity
Storage	Store dry and dark at -20°C or for long term storage at -80°C.		

\*Patents Pending