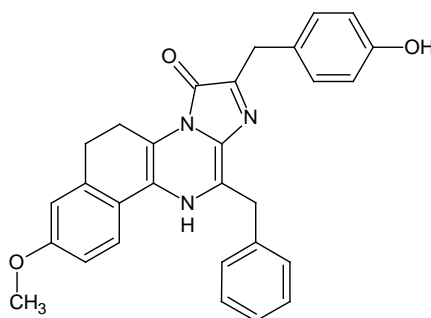


Certificate of Analysis

CAT# 369 “Prolume Purple” Methoxy-eCoelenterazine

Date of Test: 06 December 2019
 Compound ID: Methoxy-eCoelenterazine
 CAS#: Not Listed, patent pending
 IUPAC Name: Methoxy-eCoelenterazine or “Prolume Purple”
 Batch Number: L0612 (86-081-1C)
 Manufacturer: Prolume Ltd
 Molecular Formula: C₂₉H₂₅N₃O₃
 MW: 463.53 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Yellow-Orange Solid	Conforms	Visual
Purity	>95%	96.5%	HPLC
Identification	NMR is consistent with structure	Conforms	NMR
	Mass Spec is consistent with structure	Conforms	MS
Storage	Store dry and dark at -20°C or for long term storage at -80°C.		