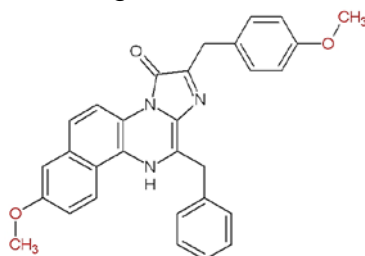


Certificate of Analysis

CAT #361 “Prolume Purple V” Methoxy-v-Coelenterazine

Date of Test: 09 June 2019
 Compound ID: Methoxy-v-Coelenterazine-Methoxy (97.7%)
 CAS#: Not Listed, patent pending
 IUPAC Name: Me-O-vCTZ-O-Me or “Prolume Purple V”
 Batch Number: L2713 (11606-038-MeO)
 Manufacturer: Prolume Ltd
 Molecular Formula: C₃₀H₂₅N₃O₃
 MW: 475.19 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Pale Yellow Solid	Conforms	Visual
Purity	>95%	97.7%	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.		