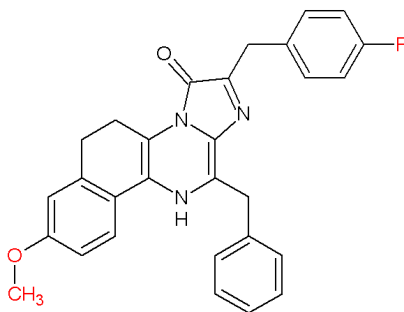


Certificate of Analysis

CAT #365 “Prolume Purple III” Methoxy-eCoelenterazineF

Date of Test: 22 May 2019
 Compound ID: Methoxy-eCoelenterazineF
 CAS#: Not Listed, patent pending
 IUPAC Name: Me-O-eCTZF or “Prolume Purple III”
 Batch Number: K1612 (05086-057)
 Manufacturer: Prolume Ltd
 Molecular Formula: C₂₉H₂₄FN₃O₂
 MW: 465.5 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Orange-Yellow Solid	Conforms	Visual
Purity	>95%	96.3%	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.		