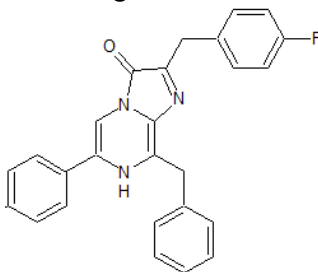


Certificate of Analysis

CAT# 348 6h-F-Coelenterazine

Date of Test: June 24, 2026
 Compound ID: 6h-F-Coelenterazine, C6d-f-coelenterazine
 CAS#: N/A
 IUPAC Name: 8-benzyl-2-(4-fluorobenzyl)-6-benzylimidazo[1,2-a]pyrazin-3(7H)-one
 Batch#: ET84801-22
 Molecular Formula: $C_{26}H_{20}FN_3O$
 Manufacturer: Prolume Ltd, patented US 9,567,623
 MW: 409.46 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Yellow Solid	Conforms	Visual
Purity	>99%	99.2%	HPLC
Identification	Spectrum is consistent with structure	Conforms	By NMR
	Spectrum is consistent with structure	Conforms	By HPLC
Storage	Store dry and dark at -20°C or for long term storage at -80°C.		