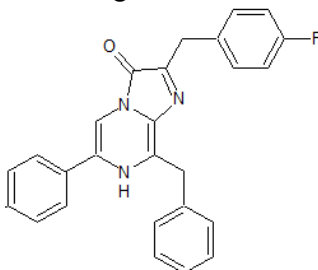


Certificate of Analysis

CAT# 3481 Water-soluble 6h-F-Coelenterazine

Date of Test: June 22, 2026
 Compound ID: 6h-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: C₂₆H₂₀FN₃O
 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. | | |