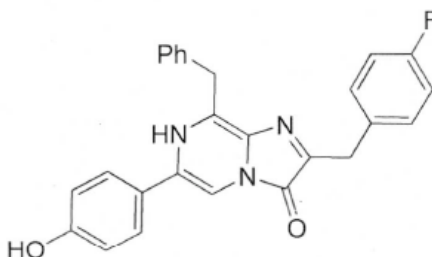


Certificate of Analysis

CAT# 345 Coelenterazine F

Date of Test: 6 September 2019
 Compound ID: Coelenterazine-Fluoride or Coelenterazine-F
 CAS#: 123437-16-1
 IUPAC Name: 8-benzyl-2-(4-fluorobenzyl)-6-(4-hydroxyphenyl)imidazo[1,2-a]pyrazin-3(7H)-one
 Batch Number: 04229-019
 Manufacturer: Prolume Ltd
 Molecular Formula: C₂₆H₂₀FN₃O₂
 MW: 425.5 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Orange-Brown Solid	Conforms	Visual
Purity	>98%	98.2%	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.		