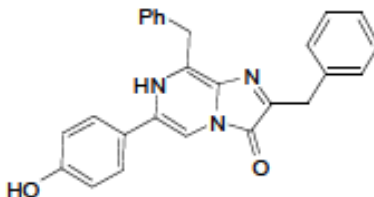


Certificate of Analysis

CAT# 301-INJ h-Coelenterazine Inject-a-Lume

Date of Test: July 22, 2019
 Compound ID: h-Coelenterazine or 2-(4-dehydroxy) Coelenterazine
 CAS#: 50909-86-9
 IUPAC Name: 2,8-dibenzyl-6-(4-hydroxyphenyl)imidazo[1,2-a]pyrazin-3(7H)-one
 Batch Number: 03611-066
 Manufacturer: Prolume Ltd
 Molecular Formula: C₂₆H₂₁N₃O₂
 MW: 407.16 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Yellow Solid	Conforms	Visual
Purity	>99%	99.7%	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.		