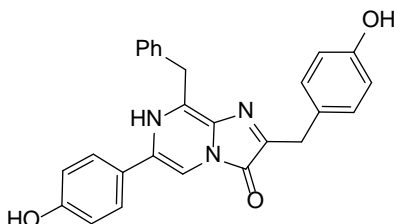


Certificate of Analysis

CAT# 3031 Water-soluble Coelenterazine native

Date of Test: June 7, 2019
 Compound ID: Coelenterazine (Native form designed for *in vivo* use)
 CAS#: 55779-48-1
 IUPAC Name: 8-benzyl-6-(4-hydroxyphenyl)-2-[(4-hydroxyphenyl)methyl]-7H-imidazo[1,2-a]pyrazin-3-one
 Batch#: 03611-052
 Molecular Formula: C₂₆H₂₁N₃O₃
 Manufacturer: Prolume Ltd
 MW: 423.46 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.		