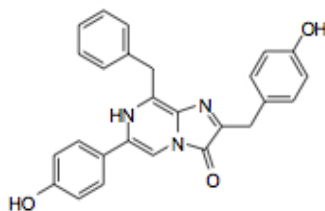


## Certificate of Analysis

### CAT# 303 Coelenterazine native

Date of Test: June 7, 2019  
 Compound ID: Coelenterazine native  
 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: C<sub>26</sub>H<sub>21</sub>N<sub>3</sub>O<sub>3</sub>  
 Manufacturer: Prolume Ltd  
 MW: 423.46 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Yellow Solid	Conforms	Visual
Purity	>97%	97.7%	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.		