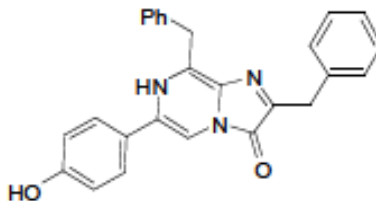


## Certificate of Analysis

### CAT# 301 h-Coelenterazine

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: C<sub>26</sub>H<sub>21</sub>N<sub>3</sub>O<sub>2</sub>  
 MW: 407.16 g/mol

Structure:



Test Items	Specification	Results	Test method
<b>Appearance</b>	Yellow Solid	Conforms	Visual
<b>Purity</b>	>99%	99.7%	HPLC
<b>Identification</b>	Spectrum is consistent with structure	Conforms	By NMR
	Spectrum is consistent with structure	Conforms	By HPLC
<b>Storage</b>	Store dry and dark at -20°C or for long term storage at -80°C.		