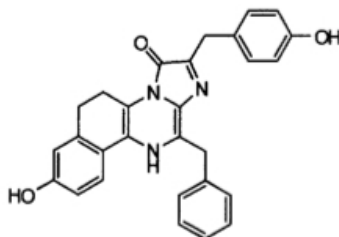


## Certificate of Analysis

### CAT# 355 e-Coelenterazine

Date of Test: 28 March 2019  
 Compound ID: e-Coelenterazine or Coelenterazine-E, eCTZ  
 CAS#: 114496-02-5  
 IUPAC Name: Benz[f]imidazo[1,2-a]quinoxalin-3(6H)-one,5,11-dihydro-8-hydroxy-2-[(4-hydroxyphenyl)methyl]-12-(phenylmethyl)  
 Batch Number: 05086-016  
 Manufacturer: Prolume Ltd  
 Molecular Formula: C<sub>28</sub>H<sub>23</sub>N<sub>2</sub>O<sub>3</sub>  
 MW: 449.5 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Yellow-Orange Solid	Conforms	Visual
Purity	>95%	95.8%	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.		