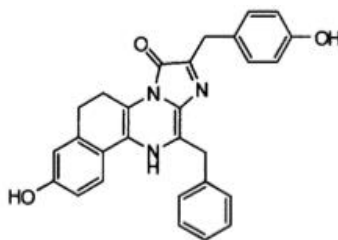


Certificate of Analysis

CAT# 3551 e-Coelenterazine-SOL

Date of Test:	24 June 2020
Compound ID:	e-Coelenterazine-SOL or water soluble Coelenterazine-E, eCTZ-SOL
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:	F2024
Manufacturer:	Prolume Ltd
Molecular Formula:	C ₂₈ H ₂₃ N ₂ O ₃
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.		