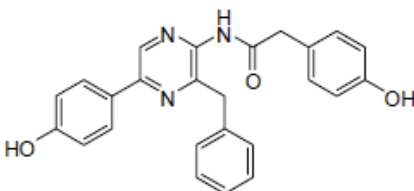


Certificate of Analysis

CAT# 350 Coelenteramide

Date of Test: 7 July 2019
 Compound ID: Coelenteramide
 CAS#: 50611-86-4
 IUPAC Name: N-[2-benzyl-6-(4-oxocyclohexa-2,5-dien-1-ylidene)-1H-pyrazin-3-yl]-2-(4hydroxyphenyl) acetamide
 Batch Number: 03647-040
 Manufacturer: Prolume Ltd
 Molecular Formula: C₂₅H₂₁N₃O₃
 MW: 411.45 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Yellow Solid	Conforms	Visual
Purity	>98%	98.47%	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.		