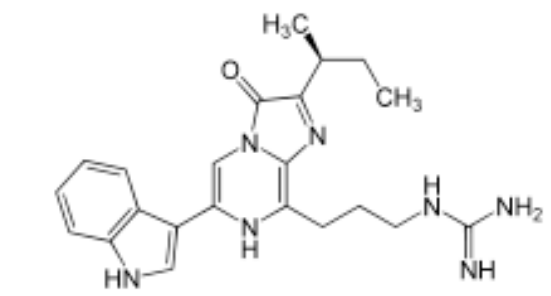


Certificate of Analysis

CAT# 305 Vargulin or Cypridina Luciferin

Date of Test: 18 December 2024
 Compound ID: Vargulin or Cypridina Luciferin
 CAS#: 7273-34-9
 IUPAC Name: 2-[3-[2-[(2S)-butan-2-yl]-6-(1H-indol-3-yl)-3-oxo-7H-imidazo[2,1-c]pyrazin-8-yl]propyl]guanidine
 Batch Number: 03784-076-H
 Manufacturer: Prolume Ltd.
 Molecular Formula: C₂₂H₂₇N₇O
 MW: 405.5 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Red to Brown Solid	Conforms	Visual
Purity	>99%	99.07%	HPLC
	Chiral purity	70% S-form	chiral-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.		