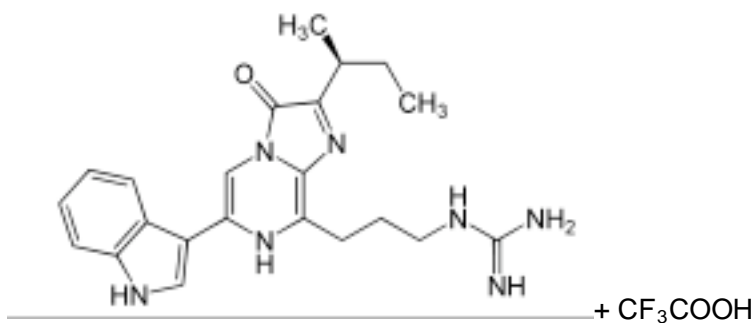


Certificate of Analysis

CAT# 305 Vargulin or Cypridina Luciferin

Date of Test: 29 June 2019
 Compound ID: Vargulin or Cypridina Luciferin - TFA salt
 CAS#: 7273-34-9
 IUPAC Name: 2-[3-[2-[(2*S*)-butan-2-yl]-6-(1*H*-indol-3-yl)-3-oxo-7*H*-imidazo[2,1-*c*]pyrazin-8-yl]propyl]guanidine trifluoroacetate
 Batch Number: 03784-076
 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
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.		