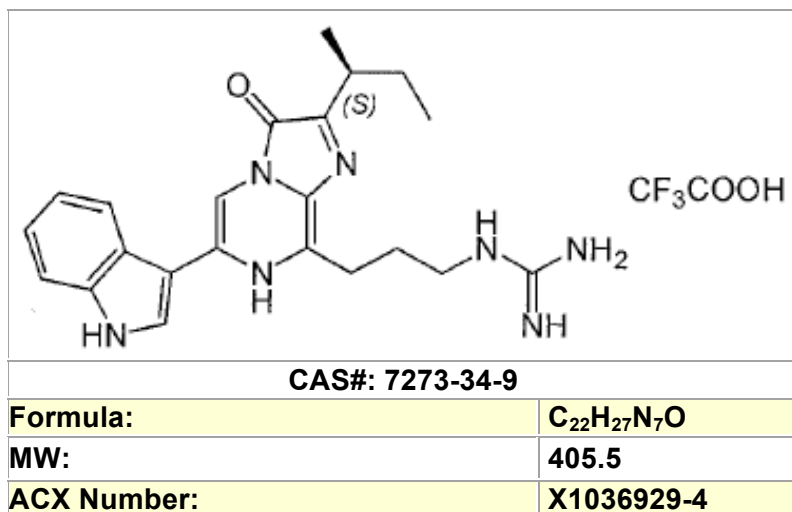


## Cypridina Luciferin (also known as Vargulin or Cypridinin)



Cypridina luciferin TFA is hundred times more unstable than native Coelenterazine.

For best results with any luciferin, dissolve immediately before use.

No matter how carefully stored, solutions will NOT KEEP without a significant loss of activity, or loss of activity that is detectable by luminometer.

To dissolve Vargulin, we suggest using Butanol, and add a trace of HCL to increase stability.

To make a short term primary storage solution: add a few microliters of 3N HCl to 1-2 ml of Butanol (or Methanol) solution until the yellow color vanishes. Then store this solution in a full stoppered tubes or under Nitrogen, Carbon Dioxide, Argon. Once dissolved if stored below -20C or -70C it may last one or a few weeks at most.

It is far more stable in n-Butanol than in Methanol, you may directly evacuate Butanol solutions under vacuum (you can't do that with Methanolic solutions).

Regardless of storage method chosen, either Vacuum or Argon, you must remove ALL the air completely; even a small amount of leftover air is as harmful as 100% air.

It is best to use Argon gas to protect the luciferin from any oxidation, keeping tight vacuum seal at -20C is often difficult (this will depend on the glassware seals that you use).

Store the dry material at -70 or in liquid Nitrogen. It has a multiyear storage life if stored as the dried TFA or 2HBr salt at low temperatures.