

## Coelenterazine h

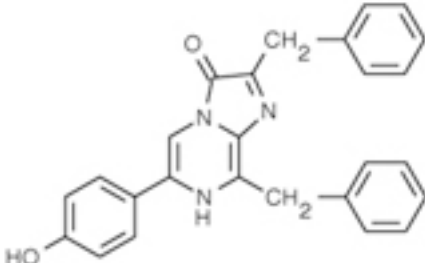
Catalog Number	Packaging Size
C239	250 µg

**Storage upon receipt:** -20°C, protected from light

### Introduction

**Coelenterazine h** is a synthetic analog of native coelenterazine and used as a bioluminescence probe for measuring  $\text{Ca}^{2+}$  concentration in cells. Coelenterazine h can generate 10 times higher luminescence intensity than native coelenterazine.

### Specifications

<b>Detection Method:</b>	Bioluminescent	
<b>Molecular Formula:</b>	$\text{C}_{26}\text{H}_{21}\text{N}_3\text{O}_2$	
<b>Molecular Weight:</b>	407.47	
<b>CAS Number:</b>	50909-86-9	
<b>Product Size:</b>	250 µg	
<b>Storage Conditions:</b>	-20°C, protected from light	
<b>Shipping Condition:</b>	Room Temperature	

### Applications

Calcium indicator

### References:

Expression of apo-aequorin during embryonic development; how much is needed for calcium imaging?  
Créton R, Steele ME, Jaffe LF  
Cell Calcium (1997) 22:439-446

Imaging Calcium Dynamics in Living Plants Cells and Tissues.  
Read ND, et al.  
Cell Biol Int (1993) 17:111-111

Chemiluminescence of Coelenterazine Analogues -Structures of Emitting Species-.  
Hirano T, et al.  
Tetrahedron Lett (1992) 39:5771-5771