

Coelenterazine

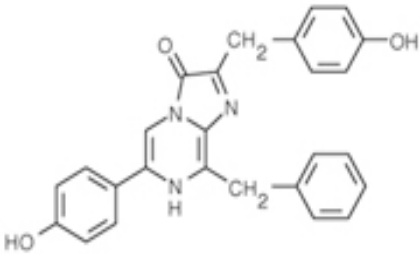
Catalog Number	Packaging Size
C238	1 mg

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

Introduction

Coelenterazine is a bioluminescence probe for measuring Ca^{2+} concentration in cells with a high sensitivity and large dynamic range. Coelenterazine (native form) is a luminescent enzyme substrate for apoaequorin and Renilla luciferase. The aequorin complex has been used as an excellent calcium indicator that can measure a broad concentration range of calcium from ~0.1 mM to >100 mM. Luminophore of the aequorin complex which is oxidized by oxygen to illuminate at 465 nm when Ca^{2+} binds to the complex.

Specifications

Detection Method:	Bioluminescent	
Molecular Formula:	$\text{C}_{26}\text{H}_{21}\text{N}_3\text{O}_3$	
Molecular Weight:	423.47	
CAS Number:	55779-48-1	
Product Size:	1 mg	
Storage Conditions:	-20°C, protected from light	
Shipping Condition:	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

A cell-based cGMP assay useful for ultra-high-throughput screening and identification of modulators of the nitric oxide/cGMP pathway.

Wunder F, Stasch JP, Hütter J, Alonso-Alija C, Hüser J, Lohrmann E