

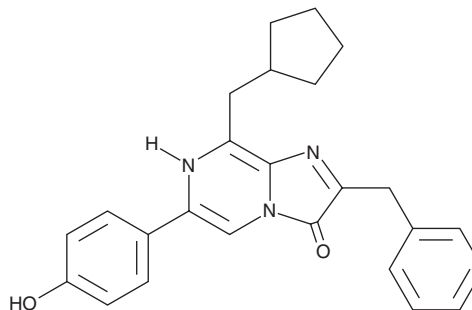
PRODUCT INFORMATION



Coelenterazine hcp

Item No. 34724

CAS Registry No.: 123437-32-1
Formal Name: 8-(cyclopentylmethyl)-6-(4-hydroxyphenyl)-2-(phenylmethyl)-imidazo[1,2-a]pyrazin-3(7H)-one
MF: C₂₅H₂₅N₃O₂
FW: 399.5
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Em. Max: 445 nm



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Coelenterazine hcp is supplied as a solid. A stock solution may be made by dissolving the coelenterazine hcp in the solvent of choice, which should be purged with an inert gas. Coelenterazine hcp is soluble in ethanol.

Description

Coelenterazine hcp is a synthetic bioluminescent luciferin that displays an emission maximum of 445 nm.¹ It has been used as a calcium indicator and substrate to quantify *Renilla* luciferase activity.¹⁻³

References

1. Sabnis, R.W. *Handbook of biological dyes and stains: Synthesis and industrial applications*. 1st ed., John Wiley & Sons, Inc., Madison, NJ, USA (2010).
2. Shimomura, O., Kishi, Y., and Inouye, S. The relative rate of aequorin regeneration from apoaequorin and coelenterazine analogues. *Biochem. J.* **296(Pt 3)**, 549-551 (1993).
3. Pichler, A., Prior, J.L., and Piwnicka-Worms, D. Imaging reversal of multidrug resistance in living mice with bioluminescence: MDR1 P-glycoprotein transports coelenterazine. *Proc. Natl. Acad. Sci. USA* **101(6)**, 1702-1707 (2004).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/22/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM