

2-(4-Dehydroxy)coelenterazine

Chemiluminescence probe of the superoxide anion and peroxynitrite

Derivative of coelenterazine (Prod. No. [ENZ-52054](#)) with ~20 times higher luminescence intensity.

Citations: 6

[View Online »](#)

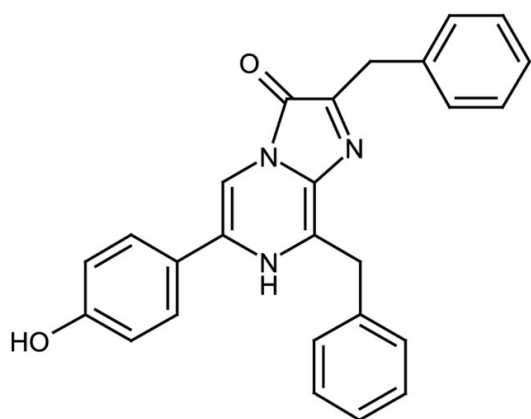
Ordering Information

[Order Online »](#)

ALX-620-062-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Keep under inert gas. Keep calcium free when stored in solution.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Coelenterazine h, Luciferin / Renilla
Appearance	Yellow to orange solid.
CAS	50909-86-9
Formula	$C_{26}H_{21}N_3O_2$
MW	407.5
Purity	≥95%
Solubility	Soluble in 100% ethanol or methanol; slightly soluble in water. Avoid DMSO as product is unstable in this solvent.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)