

# Coelenterazine (ultra pure)

## Calcium dye

Coelenterazine is useful for monitoring calcium in live cells and tissues, in reporter gene assays, for superoxide anion detection and in drug high-throughput screening applications. Coelenterazine is useful for reconstituting aequorin in cells that have been transfected with apo-aequorin cDNA. The aequorin complex comprises a 22kDa apoaequorin protein, molecular oxygen and coelenterazine. In the presence of Ca(II), coelenterazine is oxidized to coelenteramide, yielding the release of carbon dioxide and blue light. Unlike fluorescent Ca(II) indicators, Ca<sup>2+</sup>-bound aequorin can be detected without illuminating the sample, thereby eliminating interference from autofluorescence. Wavelength Maxima: Excitation 429nm, Emission 466nm

Citations: 2

[View Online »](#)

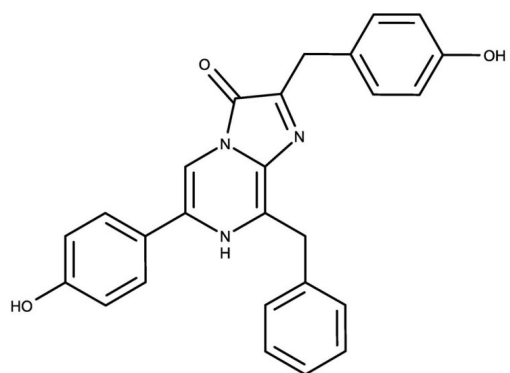
## Ordering Information

[Order Online »](#)

ENZ-52054	250µg
-----------	-------

## Manuals, SDS & CofA

[View Online »](#)



# Handling & Storage

Use/Stability	Store at -20°C, under nitrogen. Desiccation recommended. Protect material from long-term exposure to light; may be exposed to light for short periods of time. This material is unstable in DMSO, DMF, dimethylacetamide and N-methylpyrrolidone. Coelenterazine solution in methanol showed <10% decomposition when stored at room temperature in the dark for one month, and <2% decomposition when stored at -20°C in the dark, under nitrogen.
Handling	Protect from light. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	2-[(4-Hydroxyphenyl)methyl]-6-(4-hydroxyphenyl)-8-(phenylmethyl)-imidazo [1,2-a] pyrazin-3-(7H)-one
CAS	55779-48-1
Couple Type	Dye
MW	423.5
Purity	≥90% (HPLC)
Solubility	Soluble in 100% ethanol.

### Technical Info / Product Notes

This product is a member of the CELLESTIAL<sup>®</sup> product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL<sup>®</sup> reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.



ENZO LIFE SCIENCES,  
INC.  
Phone: +800.943.0420  
[usa@enzolifesciences.com](mailto: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-  
be@enzolifesciences.com](mailto:info-be@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)