

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, Ca²⁺-bound aequorin can be detected without illuminating the sample, thereby eliminating interference from autofluorescence. Wavelength Maxima: Excitation 429nm, Emission 466nm

Citations: 2

[View Online »](#)

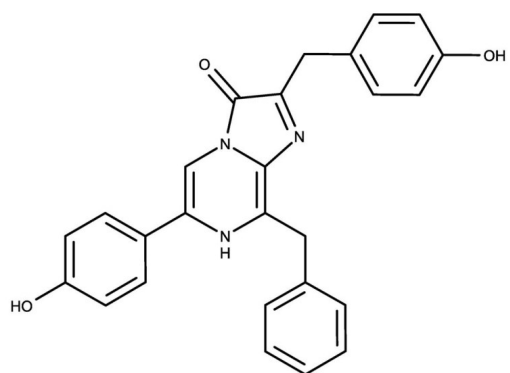
Ordering Information

[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 [®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL [®] 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: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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