

Luminol (ultra pure)

ROS dye

Luminol is a luminescent peroxidase substrate that can be used as a chemiluminescent reagent for the determination of viable mammalian cells and bacteria. Luminol also has forensic application as a presumptive test for latent blood detection and is commonly employed in detection strategies by flow injection analysis, for pharmaceutical and environmental characterization. Chemiluminescent and electrochemiluminescent probes based upon luminal are beginning to merge with biochip and microarray development as a possible substitutes for the more well established fluorescent probes. Wavelength Maxima: Excitation 355nm, Emission 411nm

Citations: 1

[View Online »](#)

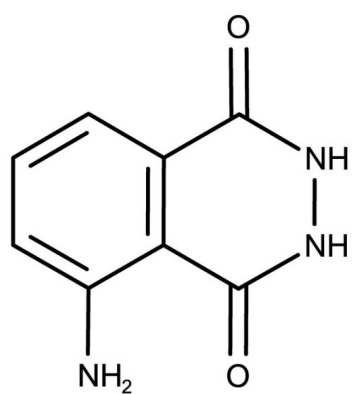
Ordering Information

[Order Online »](#)

ENZ-52354	1g
-----------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
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	3-Aminophthalhydrazide, (S)-2-(6-Hydroxy-2-benzothiazolyl)-2-thiazoline-4-carboxylic acid
CAS	521-31-3
Formula	$C_8H_7N_3O_2$
MW	177.2
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.
Technical Info / Product Notes	<p>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.</p>

Last modified: May 29, 2024



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