

Monobromobimane

(ultra pure)

Thiol dye

Monobromobimane (mBBBr) is a thiol-reactive fluorescent probe used for the determination of the redox status of low molecular weight and protein thiols in biological systems. Monobromobimane-based *in situ* derivatization results in maximal recovery of both free, reduced low molecular weight and monobromobimane-accessible protein thiols. The quantitation of the corresponding adducts of protein thiols is achieved by fluorescence spectroscopy, following protein precipitation. The reagent is particularly useful for quantifying glutathione. Wavelength Maxima: Excitation 395nm, Emission 490nm

Citations: 2

[View Online »](#)

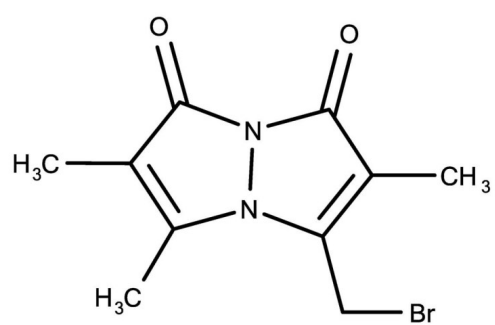
Ordering Information

[Order Online »](#)

ENZ-52501	25mg
-----------	------

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	mBBr, 3-(Bromomethyl)-2,5,6-trimethyl-1H,7H-pyrazolo(1,2- α) pyrazole-1,7-dione
CAS	71418-44-5
Formula	$C_{10}H_{11}BrN_2O_2$
MW	271.1
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](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)