

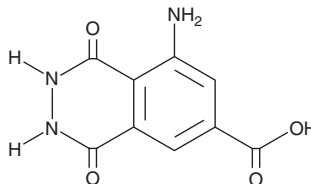
PRODUCT INFORMATION



m-carboxy Luminol

Item No. 36914

CAS Registry No.: 2221025-25-6
Formal Name: 8-amino-1,2,3,4-tetrahydro-1,4-dioxo-6-phthalazinecarboxylic acid
Synonym: *meta*-carboxy Luminol
MF: C₉H₇N₃O₄
FW: 221.2
Purity: ≥90%
Ex./Em. MAX.: 385/498.6 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

m-carboxy Luminol is supplied as a solid. A stock solution may be made by dissolving the *m*-carboxy luminol in the solvent of choice, which should be purged with an inert gas. *m*-carboxy Luminol is soluble (≥10 mg/ml) in acetonitrile:water (1:1).

Description

m-carboxy Luminol is a water-soluble chemiluminescent probe.^{1,2} It displays an emission maximum of 498.6 nm upon excitation at 385 nm.² *m*-carboxy Luminol exhibits increased chemiluminescent intensity and stability compared with luminol (Item No. 16803).^{1,2} It has been used in enzymatic activity assays and immunoassays, as well as to detect bloodstains.

References

1. Rink, S., Duerkop, A., Jacobi von Wangelin, A., *et al.* Next generation luminol derivative as powerful benchmark probe for chemiluminescence assays. *Anal. Chim. Acta* **1188**, 339161 (2021).
2. Wu, M., Liu, A., Xu, T., *et al.* Highly sensitive chemiluminescence of water-soluble luminol and its application in forensic bloodstain detection. *Anal. Methods* **17(15)**, 2934-2937 (2025).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/22/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM