

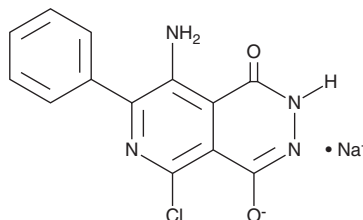
PRODUCT INFORMATION



L-012 (sodium salt)

Item No. 39908

CAS Registry No.: 143556-24-5
Formal Name: 8-amino-5-chloro-2,3-dihydro-7-phenyl-pyrido[3,4-d]pyridazine-1,4-dione, monosodium salt
MF: $C_{13}H_8ClN_4O_2 \cdot Na$
FW: 310.7
Purity: $\geq 98\%$
Em. Max: 470 nm
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

L-012 (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the L-012 (sodium salt) in the solvent of choice, which should be purged with an inert gas. L-012 (sodium salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

L-012 is a chemiluminescent probe for the detection of superoxide and a derivative of luminol (Item No. 16803).¹ It has commonly been used for the detection of NADPH oxidase-derived superoxide *in vitro* and for the detection of reactive oxygen species (ROS) and reactive nitrogen species (RNS) *in vivo*.¹⁻³ L-012 has a chemiluminescent emission maximum of 470 nm.³

References

1. Daiber, A., August, M., Baldus, S., *et al.* Measurement of NAD(P)H oxidase-derived superoxide with the luminol analogue L-012. *Free Radic. Biol. Med.* **36**(1), 101-111 (2004).
2. Kielland, A., Blom, T., Nandakumar, K.S., *et al.* *In vivo* imaging of reactive oxygen and nitrogen species in inflammation using the luminescent probe L-012. *Free Radic. Biol. Med.* **47**(6), 760-766 (2009).
3. Gao, L., Zhang, Y., Zhao, L., *et al.* An artificial metalloenzyme for catalytic cancer-specific DNA cleavage and operando imaging. *Sci. Adv.* **6**(29), eabb1421 (2020).

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, 02/21/2025

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