

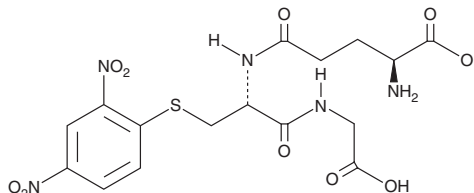
PRODUCT INFORMATION



S-(2,4-Dinitrophenyl)-Glutathione

Item No. 24660

CAS Registry No.: 26289-39-4
Formal Name: L-γ-glutamyl-S-(2,4-dinitrophenyl)-L-cysteinyl-glycine
Synonyms: Dinitrophenyl-S-Glutathione, NSC 131112, NSC 638615
MF: C₁₆H₁₉N₅O₁₀S
FW: 473.4
Purity: ≥98%
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

S-(2,4-Dinitrophenyl)-glutathione is supplied as a solid. S-(2,4-Dinitrophenyl)-glutathione is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

S-(2,4-Dinitrophenyl)-glutathione is a substrate for glutathione-S-transferase with a kinetic dissociation constant of 7 μM.¹ *In vitro*, it is transported by multidrug resistance protein 1 (MDR1) and MDR3, which are transporters that confer resistance to some chemotherapeutics.²

References

- Schramm, V.L., McCluskey, R., Emig, F.A., *et al.* Kinetic studies and active site-binding properties of glutathione S-transferase using spin-labeled glutathione, a product analogue. *J. Biol. Chem.* **259**(2), 714-722 (1984).
- Zeng, H., Liu, G., Rea, P.A., *et al.* Transport of amphipathic anions by human multidrug resistance protein 3. *Cancer Res.* **60**(17), 4779-4784 (2000).

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, 11/23/2022

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