

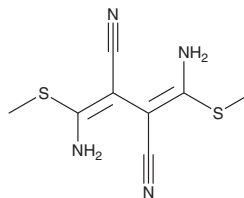
PRODUCT INFORMATION



U-0124

Item No. 34871

CAS Registry No.: 108923-79-1
Formal Name: 2,3-bis[amino(methylthio)methylene]-butanedinitrile
MF: C₈H₁₀N₄S₂
FW: 226.3
Purity: ≥95%
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

U-0124 is supplied as a solid. A stock solution may be made by dissolving the U-0124 in the solvent of choice, which should be purged with an inert gas. U-0124 is slightly soluble in DMSO and methanol.

Description

U-0124 is an inactive analog of the MEK inhibitor and AMPK activator U-0126 (Item No. 70970) that has been used as a negative control for the kinase activity of MEK.^{1,2}

References

1. Favata, M.F., Horiuchi, K.Y., Manos, E.J., *et al.* Identification of a novel inhibitor of mitogen-activated protein kinase kinase. *J. Biol. Chem.* **273**(29), 18623-18632 (1998).
2. Zipperly, M.E., Sultan, F.A., Graham, G.-E., *et al.* Regulation of dopamine-dependent transcription and cocaine action by *Gadd45b*. *Neuropsychopharmacology* **46**(4), 709-720 (2021).

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, 05/20/2024

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