

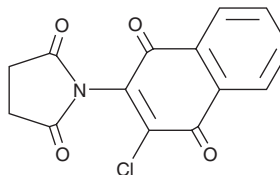
PRODUCT INFORMATION



MEK Inhibitor II

Item No. 45219

CAS Registry No.: 623163-52-0
Formal Name: 1-(3-chloro-1,4-dihydro-1,4-dioxo-2-naphthalenyl)-2,5-pyrrolidinedione
Synonyms: IMDNQ1
MF: C₁₄H₈ClNO₄
FW: 289.7
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

MEK inhibitor II is supplied as a solid. A stock solution may be made by dissolving the MEK inhibitor II in the solvent of choice, which should be purged with an inert gas. MEK inhibitor II is sparingly soluble (1-10 mg/ml) in DMSO.

Description

MEK inhibitor II is an inhibitor of MEK1 (IC₅₀ = 0.38 μM in a cell-free assay).¹ It is selective for MEK1 over ERK1 and Raf1 (IC₅₀s = 82.9 and 34.5 μM, respectively). It is also active against *T. cruzi* (IC₅₀ = 2.77 μM).²

References

1. Bakare, O., Ashendel, C.L., Peng, H., *et al.* Synthesis and MEK1 inhibitory activities of imido-substituted 2-chloro-1,4-naphthoquinones. *Bioorg. Med. Chem.* **11(14)**, 3165-3170 (2003).
2. Khraiweh, M.H., Lee, C.M., Brandy, Y., *et al.* Antitrypanosomal activities and cytotoxicity of some novel imido-substituted 1,4-naphthoquinone derivatives. *Arch. Pharm. Res.* **35(1)**, 27-33 (2012).

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, 04/29/2026

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