

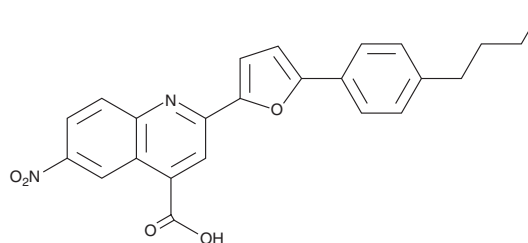
PRODUCT INFORMATION



eIF4A i28

Item No. 45394

CAS Registry No.: 2861994-92-3
Formal Name: 2-[5-(4-butylphenyl)-2-furanyl]-6-nitro-4-quinolinecarboxylic acid
Synonym: Eukaryotic Translation Initiation Factor 4A i28
MF: C₂₄H₂₀N₂O₅
FW: 416.4
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

eIF4A i28 is supplied as a solid. A stock solution may be made by dissolving the eIF4A i28 in the solvent of choice, which should be purged with an inert gas. eIF4A i28 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

eIF4A i28 is an inhibitor of eukaryotic translation initiation factor 4A (eIF4A).¹ It inhibits the ATPase and helicase activities of eIF4A with IC₅₀ values of 8.6 and 6.6 μM, respectively, in cell-free assays. eIF4A i28 reduces the viability of BJAB Burkitt lymphoma cells (EC₅₀ = 0.46 μM).

Reference

1. Zerio, C.J., Cunningham, T.A., Tulino, A.S., *et al.* Discovery of an eIF4A inhibitor with a novel mechanism of action. *J. Med. Chem.* **64**(21), 15727-15746 (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, 04/20/2026

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