

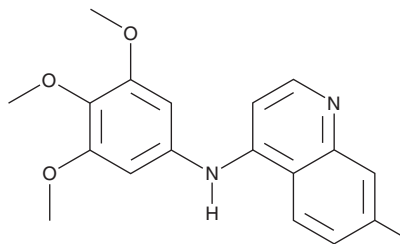
PRODUCT INFORMATION



PKN3 Inhibitor Compound 16

Item No. 41437

CAS Registry No.: 2361545-75-5
Formal Name: 7-iodo-N-(3,4,5-trimethoxyphenyl)-4-quinolinamine
Synonym: Protein Kinase N3 Inhibitor Compound 16
MF: C₁₈H₁₇IN₂O₃
FW: 436.2
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

PKN3 inhibitor compound 16 is supplied as a solid. A stock solution may be made by dissolving the PKN3 inhibitor compound 16 in the solvent of choice, which should be purged with an inert gas. PKN3 inhibitor compound 16 is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

PKN3 inhibitor compound 16 is an inhibitor of serine/threonine protein kinase N3 (IC₅₀ = 14 nM).¹ It is selective for this kinase over a panel of 234 kinases but does inhibit serine/threonine protein kinase NLK, activin receptor type 1 (ACTR1), and receptor-interacting serine/threonine kinase 2 (RIPK2) at 1 μM, as well as cyclin G-associated kinase (GAK; IC₅₀ = 74 nM in a bioluminescence resonance energy transfer (BRET) assay).^{1,2} PKN3 inhibitor compound 16 inhibits viral replication in dengue virus-infected Huh7 cells (EC₅₀ = 3.6 μM) without inducing cell death with a 50% cytotoxic concentration (CC₅₀) value of >10 μM.³

References

1. Asquith, C.R.M., Temme, L., East, M.P., *et al.* Identification of 4-anilinoquin(az)oline as a cell-active protein kinase novel 3 (PKN3) inhibitor chemotype. *ChemMedChem* **17**(12), e202200161 (2022).
2. Asquith, C.R.M., Laitinen, T., Bennett, J.M., *et al.* Design and analysis of the 4-anilinoquin(az)oline kinase inhibition profiles of GAK/SLK/STK10 using quantitative structure-activity relationships. *ChemMedChem* **15**(1), 26-49 (2020).
3. Saul, S., Pu, S.-Y., Zuercher, W.J., *et al.* Potent antiviral activity of novel multi-substituted 4-anilinoquin(az)olines. *Bioorg. Med. Chem. Lett.* **30**(16), 127284 (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, 08/30/2024

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