

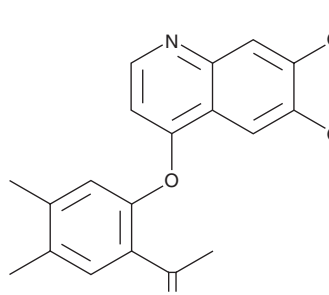
PRODUCT INFORMATION



TGF- β RI Kinase Inhibitor VII

Item No. 45225

CAS Registry No.: 666729-57-3
Formal Name: 1-[2-[(6,7-dimethoxy-4-quinolinyl)oxy]-4,5-dimethylphenyl]-ethanone
MF: $C_{21}H_{21}NO_4$
FW: 351.4
Purity: $\geq 98\%$
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

TGF- β RI kinase inhibitor VII is supplied as a solid. A stock solution may be made by dissolving the TGF- β RI kinase inhibitor VII in the solvent of choice, which should be purged with an inert gas. TGF- β RI Kinase Inhibitor VII is slightly soluble (0.1-1 mg/ml) in methanol.

Description

TGF- β RI kinase inhibitor VII is an inhibitor of activin receptor-like kinase 5 (ALK5), also known as TGF- β receptor type 1 (TGF- β R1; $IC_{50} = 0.63 \mu\text{M}$).¹ It is selective for ALK5 over c-Src ($IC_{50} = 13 \mu\text{M}$). TGF- β RI kinase inhibitor VII (0.15-5 μM) inhibits SMAD2 phosphorylation in A549 cell lysates.

Reference

1. Shimizu, T., Kimura, K., Sakai, T., *et al.* Synthesis and evaluation of [2-(4-quinolyloxy)phenyl]methanone derivatives: Novel selective inhibitors of transforming growth factor-beta kinase. *J. Med. Chem.* **51**(11), 3326-9 (2008)

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, 05/04/2026

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