

PRODUCT INFORMATION

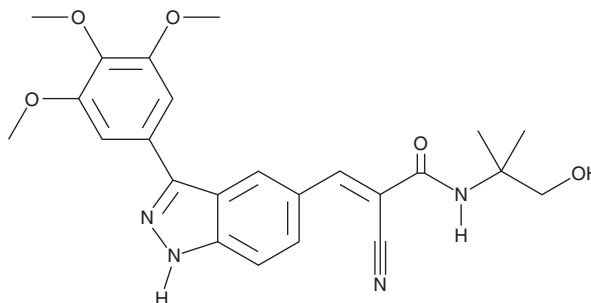


RMM 46

Item No. 35027

CAS Registry No.: 1307896-46-3
Formal Name: 2-cyano-N-(2-hydroxy-1,1-dimethylethyl)-3-[3-(3,4,5-trimethoxyphenyl)-1H-indazol-5-yl]-2-propenamide

MF: C₂₄H₂₆N₄O₅
FW: 450.5
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

RMM 46 is supplied as a solid. A stock solution may be made by dissolving the RMM 46 in the solvent of choice, which should be purged with an inert gas. RMM 46 is soluble in the organic solvent DMSO.

Description

RMM 46 is an inhibitor of p90 ribosomal S6 kinase 2 (RSK2; IC₅₀s = 12 and <2.5 nM for the wild-type and RSK2^{T493M} human enzymes, respectively).¹ It is selective for RSK2 over polo-like kinase 1 (Plk1) and NIMA-related kinase 2 (NEK2; IC₅₀s = 2,200 and 530 nM, respectively) and a panel of 26 kinases when used at a concentration of 1 μM. RMM 46 inhibits autophosphorylation of RSK2 and RSK3 induced by phorbol 12-myristate 13-acetate (PMA; Item No. 10008014) in COS-7 cells expressing the human enzymes in a concentration-dependent manner.

Reference

1. Miller, R.M., Paavilainen, V.O., Krishnan, S., *et al.* Electrophilic fragment-based design of reversible covalent kinase inhibitors. *J. Am. Chem. Soc.* **135(14)**, 5298-5301 (2013).

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, 12/09/2022

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