

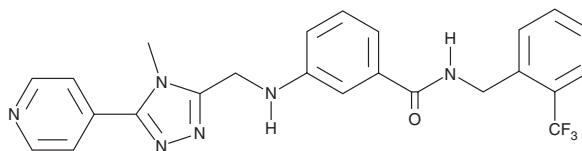
PRODUCT INFORMATION



CMPD101

Item No. 26808

CAS Registry No.: 865608-11-3
Formal Name: 3-[[[4-methyl-5-(4-pyridinyl)-4H-1,2,4-triazol-3-yl]methyl]amino]-N-[[2-(trifluoromethyl)phenyl]methyl]-benzamide
MF: C₂₄H₂₁F₃N₆O
FW: 466.5
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 247, 320 nm
Supplied as: A crystalline 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

CMPD101 is supplied as a crystalline solid. A stock solution may be made by dissolving the CMPD101 in the solvent of choice, which should be purged with an inert gas. CMPD101 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CMPD101 in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

CMPD101 is an inhibitor of G protein-coupled receptor kinase 2 (GRK2) and GRK3 (IC₅₀s = 18 and 5.4 nM, respectively).¹ It is selective for GRK2 and GRK3 over GRK1, GRK5, GRK6, and GRK7 (IC₅₀s = 3,100, 2,300, >30,000, and 25,000 nM, respectively), as well as Rho-associated kinase 2 (ROCK2) and PKCα (IC₅₀s = 1,400 and 8,100 nM, respectively). CMPD101 induces cAMP accumulation in HEK293 cells expressing human β₂-adrenergic receptors (EC₅₀ = 10 μM). In isolated human prostate strips, CMPD101 (50 μM) inhibits contractions induced by electrical field stimulation, norepinephrine, phenylephrine, endothelin-1 (Item No. 24127), and U-46619 (Item No. 16450).²

References

1. Okawa, T., Aramaki, Y., Yamamoto, M., *et al.* Design, synthesis, and evaluation of the highly selective and potent G-protein-coupled receptor kinase 2 (GRK2) inhibitor for the potential treatment of heart failure. *J. Med. Chem.* **60**(16), 6942-6990 (2017).
2. Yu, Q., Gratzke, C., Wang, Y., *et al.* Inhibition of prostatic smooth muscle contraction by the inhibitor of G protein-coupled receptor kinase 2/3, CMPD101. *Eur. J. Pharmacol.* **831**, 9-19 (2018).

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, 11/14/2022

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