

PRODUCT INFORMATION



T16A(inh)-A01

Item No. 18518

CAS Registry No.: 552309-42-9

Formal Name: 2-[(5-ethyl-1,6-dihydro-4-methyl-6-oxo-2-pyrimidinyl)thio]-N-[4-(4-methoxyphenyl)-2-thiazolyl]-acetamide

MF: $C_{19}H_{20}N_4O_3S_2$

FW: 416.5

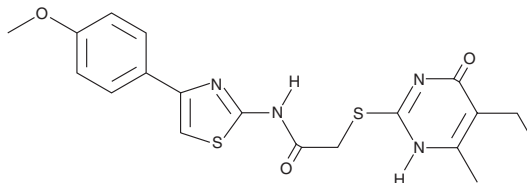
Purity: $\geq 95\%$

Supplied as: A crystalline solid

UV/Vis.: λ_{max} : 246, 274 nm

Storage: -20°C

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

T16A(inh)-A01 is supplied as a crystalline solid. A stock solution may be made by dissolving the T16A(inh)-A01 in the solvent of choice. T16A(inh)-A01 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of T16A(inh)-A01 in these solvents is approximately 5 and 10 mg/ml, respectively.

T16A(inh)-A01 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, T16A(inh)-A01 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. T16A(inh)-A01 has a solubility of approximately 0.1 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Transmembrane protein 16A (TMEM16A, also known as anoctamin-1) is a calcium-activated chloride channel (CaCC) with roles in transepithelial anion transport and smooth muscle contraction. T16A(inh)-A01 is an aminophenylthiazole that inhibits transient TMEM16A-mediated chloride currents with an IC_{50} value of $\sim 1 \mu\text{M}$.^{1,2} Its inhibitory effects are independent of voltage and do not prolong the rate of TMEM16A current deactivation.² T16A(inh)-A01 blocks CaCC activity in vascular smooth muscle cells and relaxes mouse and human blood vessels.³ It also inhibits the proliferation of pancreatic cancer and squamous carcinoma cells in culture.^{4,5}

References

1. Namkung, W., Phuan, P.W., and Verkman, A.S. TMEM16A inhibitors reveal TMEM16A as a minor component of calcium-activated chloride channel conductance in airway and intestinal epithelial cells. *J. Biol. Chem.* **286**(3), 2365-2374 (2011).
2. Bradley, E., Fedigan, S., Webb, T., et al. Pharmacological characterization of TMEM16A currents. *Channels* **8**(4), 308-320 (2014).
3. Davis, A.J., Shi, J., Pritchard, H.A.T., et al. Potent vasorelaxant activity of the TMEM16A inhibitor T16Ainh-A01. *Br. J. Pharmacol.* **168**, 773-784 (2013).
4. Mazzone, A., Eisenman, S.T., Strege, P.R., et al. Inhibition of cell proliferation by a selective inhibitor of the Ca^{2+} activated Cl^- channel, Ano1. *Biochem. Biophys. Res. Commun.* **427**(2), 248-253 (2012).
5. Duvvuri, U., Shiwarski, D.J., Xiao, D., et al. TMEM16A induces MAPK and contributes directly to tumorigenesis and cancer progression. *Cancer Res.* **72**(13), 3270-3281 (2012).

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, 12/02/2022

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