

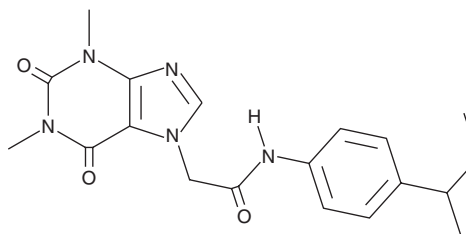
PRODUCT INFORMATION



TCS 5861528

Item No. 18380

CAS Registry No.: 332117-28-9
Formal Name: 1,2,3,6-tetrahydro-1,3-dimethyl-N-[4-(1-methylpropyl)phenyl]-2,6-dioxo-7H-purine-7-acetamide
MF: C₁₉H₂₃N₅O₃
FW: 369.4
Purity: ≥98%
UV/Vis.: λ_{max}: 248 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

TCS 5861528 is supplied as a crystalline solid. A stock solution may be made by dissolving the TCS 5861528 in the solvent of choice, which should be purged with an inert gas. TCS 5861528 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of TCS 5861528 in these solvents is approximately 0.1, 15, and 30 mg/ml, respectively.

Description

The transient receptor potential ankyrin 1 (TRPA1) ion channel integrates the nociception and transmission of diverse, potentially damaging and noxious stimuli, including cold, electrophilic compounds, divalent cations, and mechanical stimulation. TCS 5861528 is a selective TRPA1 blocker that antagonizes TRPA1-mediated calcium influx induced by allyl isothiocyanate and 4-hydroxy nonenal (Item No. 32100) with IC₅₀ values of 14.3 and 18.7 μM, respectively.^{1,2} At 30 mg/kg twice daily in rats, TCS 5861528 has been used to attenuate pain-related behavior in models of diabetic hypersensitivity and postoperative pain.^{1,3}

References

1. Wei, H., Hämäläinen, M.M., Saarnilehto, M., *et al.* Attenuation of the mechanical hypersensitivity by an antagonist of the TRPA1 ion channel in diabetic animals. *Anesthesiology* **111**, 147-154 (2009).
2. Fanger, C.M., del Camino, D., and Moran, M.M. TRPA1 as an analgesic target. *Open Drug Discov. J.* **2**, 64-70 (2010).
3. Wei, H., Karimaa, M., Korjamo, T., *et al.* Transient receptor potential ankyrin 1 ion channel contributes to guarding pain and mechanical hypersensitivity in a rat model of postoperative pain. *Anesthesiology* **117**(1), 137-148 (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, 12/14/2022

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