

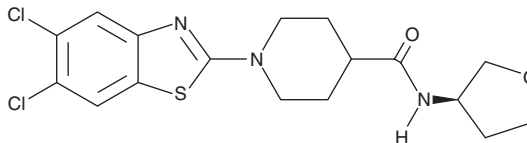
PRODUCT INFORMATION



UT-11

Item No. 38580

Formal Name: (R)-1-(5,6-dichlorobenzo[d]thiazol-2-yl)-N-(tetrahydrofuran-3-yl)piperidine-4-carboxamide
MF: C₁₇H₁₉Cl₂N₃O₂S
FW: 400.3
Purity: ≥98%
UV/Vis.: λ_{max}: 237, 284 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



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

Laboratory Procedures

UT-11 is supplied as a solid. A stock solution may be made by dissolving the UT-11 in the solvent of choice, which should be purged with an inert gas. UT-11 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of UT-11 in ethanol is approximately 3 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of UT-11 can be prepared by directly dissolving the solid in aqueous buffers. UT-11 is slightly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

UT-11 is an inhibitor of microsomal prostaglandin E synthase-1 (mPGES-1).¹ It inhibits PGE₂ production in both SK-N-AS human neuroblastoma and BV-2 mouse microglial cells (IC₅₀s = 0.1 and 2 μM, respectively). UT-11 (10 mg/kg, i.p.) reduces LPS-induced increases in a variety of cytokines, including chemokine (C-C motif) ligand 2 (CCL2), IL-6, and TNF-α, in the mouse hippocampus, but not kidney, without affecting the mRNA expression of Cox2.

Reference

- Sluter, M.N., Bhuniya, R., Yuan, X., *et al.* Novel, brain-permeable, cross-species benzothiazole inhibitors of microsomal prostaglandin E synthase-1 (mPGES-1) dampen neuroinflammation *in vitro* and *in vivo*. *ACS Pharmacol. Transl. Sci.* **6**(4), 587-599 (2023).

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, 07/12/2023

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