

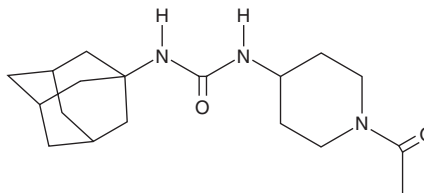
PRODUCT INFORMATION



AR-9281

Item No. 38346

CAS Registry No.: 913548-29-5
Formal Name: N-(1-acetyl-4-piperidiny)-N'-tricyclo[3.3.1.1^{3,7}]dec-1-yl-urea
Synonyms: APAU, UC1153
MF: C₁₈H₂₉N₃O₂
FW: 319.4
Purity: ≥95%
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

AR-9281 is supplied as a solid. A stock solution may be made by dissolving the AR-9281 in the solvent of choice, which should be purged with an inert gas. AR-9281 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AR-9281 in ethanol is approximately 30 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

AR-9281 is an inhibitor of soluble epoxide hydrolase (sEH; IC₅₀ = 7 nM for the human enzyme).¹ It reduces LPS-induced increases in plasma levels of dihydroxyeicosatrienoic acids (DiHETs), prostaglandin E₂ (PGE₂; Item No. 14010), and PGD₂ (Item No. 12010) in rats when administered at a dose of 3 mg/kg.² AR-9281 (1 and 3 mg/kg) inhibits LPS-induced allodynia in a rat model of inflammatory pain and increases the paw withdrawal threshold in a rat model of diabetic neuropathic pain induced by streptozotocin (Item No. 13104).

References

1. Jones, P.D., Tsai, H.J., Do, Z.N., *et al.* Synthesis and SAR of conformationally restricted inhibitors of soluble epoxide hydrolase. *Bioorg. Med. Chem. Lett.* **16(19)**, 5212-5216 (2006).
2. Wagner, K., Inceoglu, B., Dong, H., *et al.* Comparative efficacy of 3 soluble epoxide hydrolase inhibitors in rat neuropathic and inflammatory pain models. *Eur. J. Pharmacol.* **700(1-3)**, 93-101 (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, 08/14/2023

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