

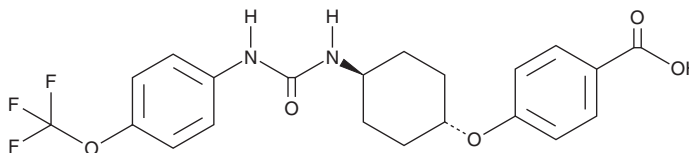
PRODUCT INFORMATION



t-TUCB

Item No. 36384

CAS Registry No.: 948304-40-3
Formal Name: 4-[[*trans*-4-[[[4-(trifluoromethoxy)phenyl]amino]carbonyl]amino]cyclohexyl]oxy]-benzoic acid
MF: $C_{21}H_{21}F_3N_2O_5$
FW: 438.4
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 247 nm
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

t-TUCB is supplied as a solid. A stock solution may be made by dissolving the t-TUCB in the solvent of choice, which should be purged with an inert gas. t-TUCB is soluble in the organic solvent DMSO.

Description

t-TUCB is a soluble epoxide hydrolase (sEH) inhibitor (IC_{50} s = 0.4, 3.1, and 11.9 nM for the human, mouse, and rat enzymes, respectively).¹ It also inhibits fatty acid amide hydrolase (FAAH; IC_{50} = 260 nM for the human enzyme). t-TUCB (3 mg/kg) increases the mechanical paw withdrawal threshold in a mouse model of diabetic neuropathic pain induced by streptozotocin (Item No. 13104).¹ It prevents EKG abnormalities induced by the β_1 - and β_2 -adrenergic receptor agonist isoproterenol (Item No. 15592), as well as reduces isoproterenol-induced increases in infarct size and heart weight, in a rat model of myocardial ischemia when administered at doses of 3, 10, and 30 mg/kg.²

References

1. Sasso, O., Wagner, K., Morisseau, C., *et al.* Peripheral FAAH and soluble epoxide hydrolase inhibitors are synergistically antinociceptive. *Pharmacol. Res.* **97**, 7-15 (2015).
2. Ayush, S., Krishnamurthy, P.T., Thomas, P., *et al.* Soluble epoxide hydrolase inhibitor, t-TUCB, protects against myocardial ischaemic injury in rats. *J. Pharm. Pharmacol.* **66(9)**, 1251-1258 (2014).

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, 05/08/2023

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