

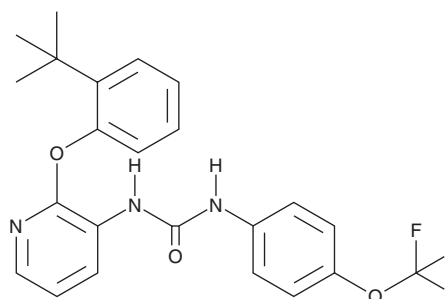
PRODUCT INFORMATION



BPTU

Item No. 29444

CAS Registry No.: 870544-59-5
Formal Name: N-[2-[2-(1,1-dimethylethyl)phenoxy]-3-pyridinyl]-N'-[4-(trifluoromethoxy)phenyl]-urea
MF: C₂₃H₂₂F₃N₃O₃
FW: 445.4
Purity: ≥98%
UV/Vis.: λ_{max}: 260, 288 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

BPTU is supplied as a solid. A stock solution may be made by dissolving the BPTU in the solvent of choice, which should be purged with an inert gas. BPTU is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BPTU in ethanol and DMF is approximately 15 mg/ml and approximately 25 mg/ml in DMSO.

BPTU is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BPTU should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BPTU has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BPTU is an allosteric antagonist of the purinergic P2Y₁ receptor (K_i = 6 nM for the human receptor).^{1,2} BPTU is selective for P2Y₁ over P2Y₂, P2Y₆, P2Y₁₁, P2Y₁₂, and P2Y₁₄ receptors (K_is = ≥3,500 nM).¹ It inhibits platelet aggregation in human platelet-rich plasma (hPRP; IC₅₀ = 2.1 μM). BPTU reduces thrombus weight by 68% in a rat model of iron chloride-induced arterial thrombosis and prolongs provoked cuticle and mesenteric bleeding time in rats when administered intravenously as a 10 mg/kg bolus dose followed by a 10 mg/kg per hour infusion. It also inhibits contractions induced by electrical field stimulation in isolated muscle strips from rat and mouse colon (EC₅₀s = 0.5 and 0.1 μM, respectively) and antrum (EC₅₀s = 9 and 0.3 μM, respectively).²

References

1. Chao, H., Turdi, H., Herpin, T.F., *et al.* Discovery of 2-(phenoxy-pyridine)-3-phenylureas as small molecule P2Y₁ antagonists. *J. Med. Chem.* **56**(4), 1704-1714 (2013).
2. Mañé, N., Jiménez-Sábado, V., and Jiménez, M. BPTU, an allosteric antagonist of P2Y₁ receptor, blocks nerve mediated inhibitory neuromuscular responses in the gastrointestinal tract of rodents. *Neuropharmacology* **110**(Pt A), 376-385 (2016).

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, 11/07/2022

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