

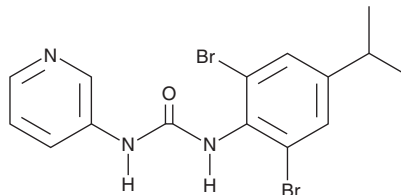
PRODUCT INFORMATION



BX-430

Item No. 32868

CAS Registry No.: 688309-70-8
Formal Name: N-[2,6-dibromo-4-(1-methylethyl)phenyl]-N'-3-pyridinyl-urea
Synonym: Blocker of P2X₄ Compound 3.0
MF: C₁₅H₁₅Br₂N₃O
FW: 413.1
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Light sensitive



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

Laboratory Procedures

BX-430 is supplied as a solid. A stock solution may be made by dissolving the BX-430 in the solvent of choice, which should be purged with an inert gas. BX-430 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

BX-430 is a noncompetitive allosteric antagonist of the purinergic P2X₄ receptor (IC₅₀s = 0.54 and 1.89 μM for the human and *Xenopus* receptors, respectively).¹ BX-430 is selective for the human P2X₄ receptor over the human P2X₁, P2X₂, P2X₅, P2X₃, and P2X₇ receptors at 5 μM and over mouse and rat P2X₄ receptors at 10 μM.

Reference

1. Ase, A.R., Honson, N.S., Zaghdane, H., *et al.* Identification and characterization of a selective allosteric antagonist of human P2X₄ receptor channels. *Mol. Pharmacol.* **87**(4), 606-616 (2015).

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, 10/06/2022

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