

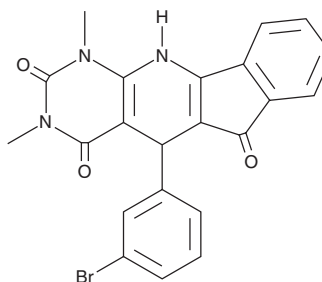
PRODUCT INFORMATION



BPIPP

Item No. 21675

CAS Registry No.: 325746-94-9
Formal Name: 5-(3-bromophenyl)-5,11-dihydro-1,3-dimethyl-1H-indeno[2',1':5,6]pyrido[2,3-d]pyrimidine-2,4,6(3H)-trione
MF: C₂₂H₁₆BrN₃O₃
FW: 450.3
Purity: ≥98%
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

BPIPP is supplied as a solid. A stock solution may be made by dissolving the BPIPP in the solvent of choice, which should be purged with an inert gas. BPIPP is slightly soluble in DMSO and methanol.

Description

BPIPP is a non-competitive inhibitor of guanylyl and adenylyl cyclases that suppresses cGMP accumulation induced by stable toxin (STa), guanylin, atrial natriuretic peptide (Item Nos. 24539 | 24276), and C-type natriuretic peptide (Item No. 24401) *in vitro* (IC₅₀s = 3.4-11.2 μM).¹ It inhibits chloride efflux induced by STa, forskolin (Item No. 11018), cholera toxin (Item No. 19654), and isoproterenol (Item No. 15592) in T84 cells but has no effect on basal chloride efflux. BPIPP (10 or 50 μM) suppresses STa-induced liquid secretion in a rabbit intestinal loop model of toxin-induced secretory diarrhea.

Reference

1. Kots, A.Y., Choi, B.-K., Estrella-Jimenez, M.E., *et al.* Pyridopyrimidine derivatives as inhibitors of cyclic nucleotide synthesis: Application for treatment of diarrhea. *Proc. Natl. Acad. Sci. U.S.A.* **105**(24), 8440-8445 (2008).

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, 11/22/2022

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