

PRODUCT INFORMATION

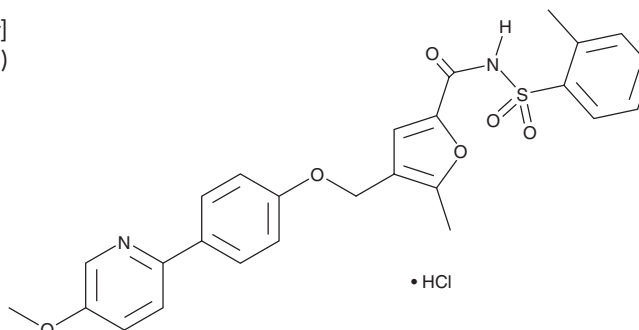


BGC 20-1531 (hydrochloride)

Item No. 19742

CAS Registry No.: 1962928-26-2
Formal Name: 4-[[4-(5-methoxy-2-pyridinyl)phenoxy]methyl]-5-methyl-N-[(2-methylphenyl)sulfonyl]-2-furancarboxamide, monohydrochloride

Synonym: PGN 1531
MF: C₂₆H₂₄N₂O₆S • HCl
FW: 529.0
Purity: ≥98%
UV/Vis.: λ_{max}: 202, 273 nm
Supplied as: A crystalline 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

BGC 20-1531 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the BGC 20-1531 (hydrochloride) in the solvent of choice. BGC 20-1531 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of BGC 20-1531 (hydrochloride) in these solvents is approximately 1 mg/ml.

Description

BGC 20-1531 is a potent and selective antagonist of the prostaglandin E₂ (PGE₂; Item No. 14010) receptor subtype 4 (EP₄) with K_i values of 11.7, >10,000, and >10,000 nM for EP₄, EP₂, and EP₃, respectively.¹ It is selective for EP₄, exhibiting <50% inhibition at 47 ion channels, cell-surface transporters, enzymes, and nuclear receptors at a concentration of 10 μM. BGC 20-1531 antagonizes PGE₂-induced cAMP accumulation in a dose-dependent manner in HEK293 EBNA cells that stably express human EP₄ receptors. It also reverses PGE₂-induced vasorelaxation (K_b = 15.85 nM) in human middle cerebral and middle meningeal arterial rings. BGC 20-1531 prevents PGE₂-induced increases in carotid blood flow and decreases in carotid resistance with an ID₅₀ value of 5 mg/kg in dogs.

Reference

1. Maubach, K.A., Davis, R.J., Clark, D.E., *et al.* BGC20-1531, a novel, potent and selective prostanoid EP₄ receptor antagonist: A putative new treatment for migraine headache. *Br. J. Pharmacol.* **156**(2), 316-327 (2009).

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, 10/17/2022

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