

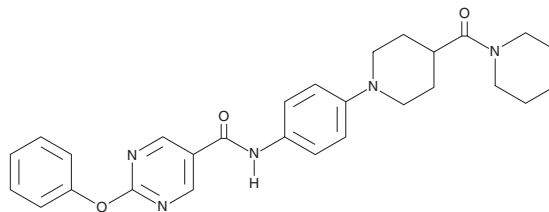
PRODUCT INFORMATION



TFC 007

Item No. 17971

CAS Registry No.: 927878-49-7
Formal Name: N-[4-[4-(4-morpholinylcarbonyl)-1-piperidinyl]phenyl]-2-phenoxy-5-pyrimidinecarboxamide
MF: C₂₇H₂₉N₅O₄
FW: 487.6
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 311 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

TFC 007 is supplied as a crystalline solid. A stock solution may be made by dissolving the TFC 007 in the solvent of choice, which should be purged with an inert gas. TFC 007 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of TFC 007 in these solvents is approximately 1 and 10 mg/ml, respectively.

Description

TFC 007 is an inhibitor of hematopoietic prostaglandin D synthase (H-PGDS; IC₅₀ = 0.083 μM).¹ It is selective for H-PGDS over lipocalin PGDS (L-PGDS) and microsomal PGE₂ synthase (mPGES) at 100 μM. It is also selective for H-PGDS over COX-1, COX-2, 5-lipoxygenase (5-LO), leukotriene C₄ (LTC₄) synthase, and thromboxane (TX) synthase, as well as the histamine H₁ and cysteinyl leukotriene 1 (CysLT₁) receptors, at 10 μM. TFC 007 (30 mg/kg) inhibits antigen-induced increases in the production of PGD₂ (Item No. 12010) in nasal cavity lavage fluid and reduces nasal blockage in a guinea pig model of allergic rhinitis.

Reference

1. Nabe, T., Kuriyama, Y., Mizutani, N., *et al.* Inhibition of hematopoietic prostaglandin D synthase improves allergic nasal blockage in guinea pigs. *Prostaglandins Other Lipid Mediat.* **95(1-4)**, 27-34 (2011).

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, 12/01/2021

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