

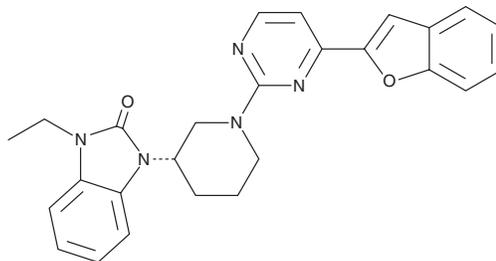
PRODUCT INFORMATION



TC-F 2

Item No. 42395

CAS Registry No.: 1304778-15-1
Formal Name: 1-[(3S)-1-[4-(2-benzofuranyl)-2-pyrimidinyl]-3-piperidinyl]-3-ethyl-1,3-dihydro-2H-benzimidazol-2-one
MF: C₂₆H₂₅N₅O₂
FW: 439.5
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

TC-F 2 is supplied as a solid. A stock solution may be made by dissolving the TC-F 2 in the solvent of choice, which should be purged with an inert gas. TC-F 2 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

TC-F 2 is an inhibitor of fatty acid amide hydrolase (FAAH; IC₅₀s = 0.028 and 0.1 μM for the human and rat enzymes, respectively).¹

Reference

1. Min, X., Thibault, S.T., Porter, A.C., et al. Discovery and molecular basis of potent noncovalent inhibitors of fatty acid amide hydrolase (FAAH). Proc. Natl. Acad. Sci. U S A 108(18), 7379-7384 (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, 12/06/2024

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