

PRODUCT INFORMATION



BAY-885

Item No. 34279

CAS Registry No.: 2307249-33-6
Formal Name: [2-amino-4-(trifluoromethoxy)phenyl][4-[7-(4-methyl-1-piperazinyl)pyrido[3,2-d]pyrimidin-4-yl]-1-piperidinyl]-methanone

MF: C₂₅H₂₈F₃N₇O₂

FW: 515.5

Purity: ≥98%

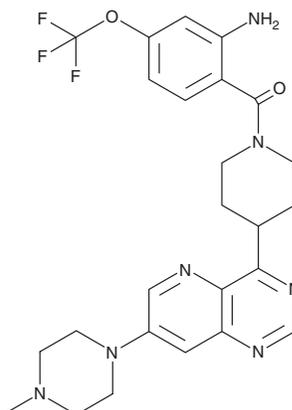
UV/Vis.: λ_{max}: 257 nm

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

BAY-885 is supplied as a solid. A stock solution may be made by dissolving the BAY-885 in the solvent of choice, which should be purged with an inert gas. BAY-885 is soluble in the organic solvent ethanol.

Description

BAY-885 is an inhibitor of ERK5 (IC₅₀ = 0.035 μM).¹ It inhibits ERK5-induced transcription in SN12C-MEF2 reporter cells (IC₅₀ = 0.12 μM).

Reference

1. Nguyen, D., Lemos, C., Wortmann, L., *et al.* Discovery and characterization of the potent and highly selective (piperidin-4-yl)pyrido[3,2-d]pyrimidine based *in vitro* probe BAY-885 for the kinase ERK5. *J. Med. Chem.* **62**(2), 928-940 (2019).

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, 11/18/2022

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