

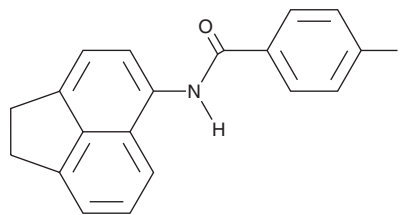
PRODUCT INFORMATION



Ebio1

Item No. 40604

CAS Registry No.: 339287-36-4
Formal Name: N-(1,2-dihydro-5-acenaphthylenyl)-4-fluoro-benzamide
MF: C₁₉H₁₄FNO
FW: 291.3
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

Ebio1 is supplied as a solid. A stock solution may be made by dissolving the ebio1 in the solvent of choice, which should be purged with an inert gas. Ebio1 is soluble in DMSO and methanol.

Description

Ebio1 is an activator of the voltage-gated potassium channel K_v7.2.¹ It enhances voltage-induced current amplitude in CHO cells expressing recombinant human K_v7.2 when used at a concentration of 10 μM.

Reference

1. Zhang, S., Ma, D., Wang, K., *et al.* A small-molecule activation mechanism that directly opens the KCNQ2 channel. *Nat. Chem. Biol.* (2024)

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, 03/06/2024

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