

PRODUCT INFORMATION



ANA-12

Item No. 14309

CAS Registry No.: 219766-25-3
Formal Name: N-[2-[[[(hexahydro-2-oxo-1H-azepin-3-yl)amino]carbonyl]phenyl]-benzo[b]thiophene-2-carboxamide

MF: C₂₂H₂₁N₃O₃S

FW: 407.5

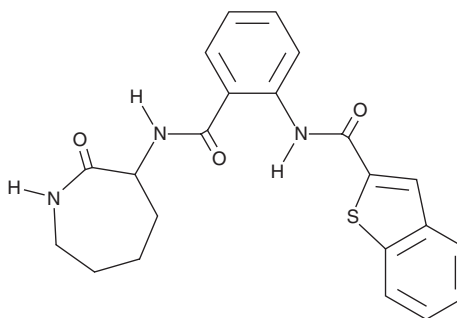
Purity: ≥98%

UV/Vis.: λ_{max}: 247, 299, 315 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

ANA-12 is supplied as a crystalline solid. A stock solution may be made by dissolving the ANA-12 in the solvent of choice, which should be purged with an inert gas. ANA-12 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ANA-12 in these solvents is approximately 5 mg/ml.

Description

ANA-12 is a TrkB receptor antagonist that prevents activation of the receptor by BDNF with IC₅₀ values of 45.6 nM and 41.1 μM for the high and low affinity sites of the receptor, respectively.¹ It does not affect TrkA or TrkC in an assay of neurite outgrowth.¹ Mice administered ANA-12 demonstrated reduced anxiety- and depression-related behaviors on a variety of tests predictive of anxiolytic and antidepressant properties in humans.¹

Reference

1. Cazorla, M., Prémont, J., Mann, A., *et al.* Identification of a low-molecular weight TrkB antagonist with anxiolytic and antidepressant activity in mice. *J. Clin. Invest.* **121**(5), 1846-1857 (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, 10/27/2022

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