

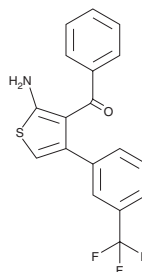
PRODUCT INFORMATION



VCP171

Item No. 30309

CAS Registry No.: 1018830-99-3
Formal Name: [2-amino-4-[3-(trifluoromethyl)phenyl]-3-thienyl]phenyl-methanone
MF: C₁₈H₁₂F₃NOS
FW: 347.4
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 241 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

VCP171 is supplied as a crystalline solid. A stock solution may be made by dissolving the VCP171 in the solvent of choice, which should be purged with an inert gas. VCP171 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of VCP171 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

VCP171 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VCP171 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. VCP171 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

VCP171 is a positive allosteric modulator of adenosine A₁ receptors (EC₅₀ = 15.8 μM in a kinetic assay measuring agonist dissociation).¹ It reduces AMPA receptor-mediated evoked excitatory postsynaptic currents (eEPSCs) in lamina I and lamina II neurons in a rat model of neuropathic pain (EC₅₀s = 1.995 and 0.251 μM, respectively) to a greater extent than in sham control animals (EC₅₀s = 2.512 and 0.631 μM, respectively).²

References

1. Aurelio, L., Figler, H., Flynn, B.L., *et al.* 5-Substituted 2-aminothiophenes as A₁ adenosine receptor allosteric enhancers. *Bioorg. Med. Chem.* **16**(3), 1319-1327 (2008).
2. Imlach, W.L., Bhola, R.F., May, L.T., *et al.* A positive allosteric modulator of the adenosine A₁ receptor selectively inhibits primary afferent synaptic transmission in a neuropathic pain model. *Mol. Pharm.* **88**(3), 460-468 (2015).

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, 08/15/2025

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