

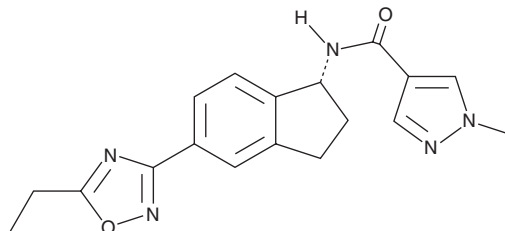
PRODUCT INFORMATION



Aficamten

Item No. 36304

CAS Registry No.: 2364554-48-1
Formal Name: N-[(1R)-5-(5-ethyl-1,2,4-oxadiazol-3-yl)-2,3-dihydro-1H-inden-1-yl]-1-methyl-1H-pyrazole-4-carboxamide
Synonyms: CK-274, CK-3773274
MF: C₁₈H₁₉N₅O₂
FW: 337.4
Purity: ≥95%
UV/Vis.: λ_{max}: 232, 248 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Aficamten is supplied as a crystalline solid. A stock solution may be made by dissolving the aficamten in the solvent of choice, which should be purged with an inert gas. Aficamten is soluble in the organic solvent DMSO.

Description

Aficamten is an inhibitor of cardiac myosin ATPase (IC₅₀ = 1.26 μM for the bovine enzyme).¹ It reduces fractional shortening in isolated rat cardiac ventricular myocytes (IC₅₀ = 7.9 μM) and in rats in a dose-dependent manner.

Reference

1. Chuang, C., Collibee, S., Ashcraft, L., *et al.* Discovery of aficamten (CK-274), a next-generation cardiac myosin inhibitor for the treatment of hypertrophic cardiomyopathy. *J. Med. Chem.* **64**(19), 14142-14152 (2021).

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, 03/04/2022

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