

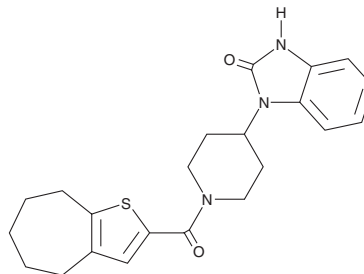
PRODUCT INFORMATION



GSK1702934A

Item No. 34576

CAS Registry No.: 924377-85-5
Formal Name: 1,3-dihydro-1-[1-[(5,6,7,8-tetrahydro-4H-cyclohepta[b]thien-2-yl)carbonyl]-4-piperidinyl]-2H-benzimidazol-2-one
MF: C₂₂H₂₅N₃O₂S
FW: 395.5
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

GSK1702934A is supplied as a solid. A stock solution may be made by dissolving the GSK1702934A in the solvent of choice, which should be purged with an inert gas. GSK1702934A is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

GSK1702934A is an activator of transient receptor potential canonical 3 (TRPC3) and TRPC6.¹ It induces TRPC3- and TRPC6-dependent currents in whole-cell patch-clamp assays using HEK293 cells expressing the human channels (EC₅₀s = 0.08 and 0.44 μM, respectively). GSK1702934A enhances contractility and induces arrhythmias in hearts isolated from TRPC3-overexpressing, but not wild-type, mice.²

References

1. Xu, X., Lozinskaya, I., Costell, M., *et al.* Characterization of small molecule TRPC3 and TRPC6 agonist and antagonists. *Biophys. J.* **104**(2), 454A (2013).
2. Doleschal, B., Primessnig, U., Wölkart, G., *et al.* TRPC3 contributes to regulation of cardiac contractility and arrhythmogenesis by dynamic interaction with NCX1. *Cardiovasc. Res.* **106**(1), 163-173 (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, 12/15/2022

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