

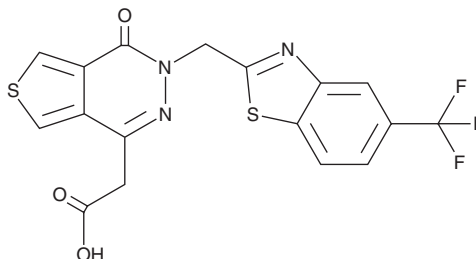
PRODUCT INFORMATION



AT-007

Item No. 44265

CAS Registry No.: 2170729-29-8
Formal Name: 3,4-dihydro-4-oxo-3-[[5-(trifluoromethyl)-2-benzothiazolyl]methyl]-thieno[3,4-d]pyridazine-1-acetic acid
Synonym: Goverestat
MF: C₁₇H₁₀F₃N₃O₃S₂
FW: 425.4
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

AT-007 is supplied as a solid. A stock solution may be made by dissolving the AT-007 in the solvent of choice, which should be purged with an inert gas. AT-007 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

AT-007 is an inhibitor of aldose reductase (IC₅₀ = 100 pM).¹ It reduces high glucose-induced senescence in isolated human keratinocytes when used at a concentration of 0.2 nM.²

References

1. Zhu, Y., Lobato, A.G., Rebelo, A.P., *et al.* Sorbitol reduction via govorestat ameliorates synaptic dysfunction and neurodegeneration in sorbitol dehydrogenase deficiency. *JCI Insight* **8**(10), e164954 (2023).
2. Yepuri, G., Kancharla, K., Perfetti, R., *et al.* The aldose reductase inhibitors AT-001, AT-003 and AT-007 attenuate human keratinocyte senescence. *Front. Aging* **5**, 1466281 (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, 08/26/2025

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