

PRODUCT INFORMATION

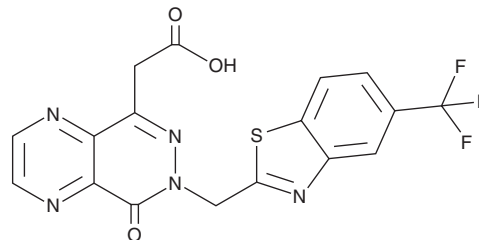


AT-001

Item No. 36105

CAS Registry No.: 1355612-71-3
Formal Name: 7,8-dihydro-8-oxo-7-[[5-(trifluoromethyl)-2-benzothiazolyl]methyl]-pyrazino[2,3-d]pyridazine-5-acetic acid

Synonym: Caficrestat
MF: C₁₇H₁₀F₃N₅O₃S
FW: 421.4
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 257 nm
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-001 is supplied as a solid. A stock solution may be made by dissolving the AT-001 in the solvent of choice, which should be purged with an inert gas. AT-001 is soluble in DMSO.

Description

AT-001 is an inhibitor of aldose reductase (IC₅₀ = 28.9 pM).¹ It reduces infarct area in streptozotocin-induced diabetic mice expressing human aldose reductase in a model of cardiomyopathy induced by ischemia-reperfusion injury when administered at a dose of 40 mg/kg.²

References

1. Wasmuth, A., Landry, D., Deng, S.X., *et al.* Aldose reductase inhibitors and uses thereof. *Trustees of Columbia University. WO2014/113380A1* (2014).
2. Perfetti, R., Rowell, P., Shendelman, S., *et al.* Preclinical and clinical proof of concept for metabolic intervention in diabetic cardiomyopathy. *J. Card. Fail.* **25(8)**, S77 (2019).

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, 12/19/2022

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