

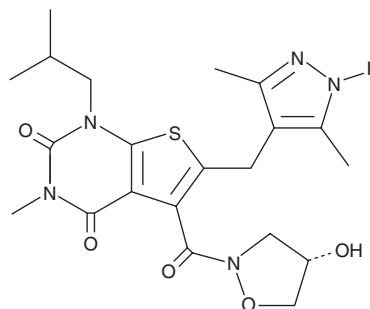
PRODUCT INFORMATION



AR-C 155858

Item No. 14884

CAS Registry No.: 496791-37-8
Formal Name: 6-[(3,5-dimethyl-1H-pyrazol-4-yl)methyl]-5-[[[(4S)-4-hydroxy-2-isoxazolidinyl]carbonyl]-3-methyl-1-(2-methylpropyl)-thieno[2,3-d]pyrimidine-2,4(1H,3H)-dione
MF: C₂₁H₂₇N₅O₅S
FW: 461.5
Purity: ≥95%
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

AR-C 155858 is supplied as a solid. A stock solution may be made by dissolving the AR-C 155858 in the solvent of choice, which should be purged with an inert gas. AR-C 155858 is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

AR-C 155858 is an inhibitor of monocarboxylate transporter 1 (MCT1) and MCT2 (K_i s = 2.3 and <10 nM, respectively, in *Xenopus* oocytes expressing the human receptors).¹ It is selective for MCT1 and MCT2 over MCT4 (K_i = >10 μM). AR-C 155858 inhibits L-lactate uptake in and proliferation of murine 4T1 breast cancer cells (IC_{50} s = 25 and 20.2 nM, respectively).²

References

- Ovens, M.J., Davies, A.J., Wilson, M.C., *et al.* AR-C155858 is a potent inhibitor of monocarboxylate transporters MCT1 and MCT2 that binds to an intracellular site involving transmembrane helices 7-10. *Biochem. J.* **425**(3), 523-530 (2010).
- Guan, X., Bryniarski, M.A., and Morris, M.E. *In vitro* and *in vivo* efficacy of the monocarboxylate transporter 1 inhibitor AR-C155858 in the murine 4T1 breast cancer tumor model. *AAPS J.* **21**(1), 3 (2018).

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, 12/01/2022

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