

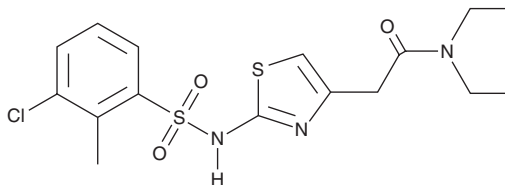
PRODUCT INFORMATION



BVT 14225

Item No. 41224

CAS Registry No.: 376638-65-2
Formal Name: 2-[[[3-chloro-2-methylphenyl)sulfonyl]amino]-N,N-diethyl-4-thiazoleacetamide
MF: C₁₆H₂₀ClN₃O₃S₂
FW: 401.9
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

BVT 14225 is supplied as a solid. A stock solution may be made by dissolving the BVT 14225 in the solvent of choice, which should be purged with an inert gas. BVT 14225 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

BVT 14225 is an inhibitor of 11 β -hydroxysteroid dehydrogenase type 1 (11 β -HSD1; IC₅₀S = 52 and 284 nM for the human and mouse enzymes, respectively).¹ It is selective for human 11 β -HSD1 over human 11 β -HSD2 (IC₅₀ = >10,000 nM). It inhibits 11 β -HSD1 by 55% in HEK293 cells expressing human *HSD11B1*, the gene encoding 11 β -HSD1, when used at a concentration of 10 μ M.² BVT 14225 (50 mg/kg) reduces blood glucose levels in a rat model of diabetes induced by streptozotocin (Item No. 13104) and nicotinamide (Item No. 11127).

References

1. Barf, T., Vallgård, J., Emond, R., *et al.* Arylsulfonamidothiazoles as a new class of potential antidiabetic drugs. Discovery of potent and selective inhibitors of the 11 β -hydroxysteroid dehydrogenase type 1. *J. Med. Chem.* **45**(18), 3813-3815 (2002).
2. Navarrete-Vázquez, G., Morales-Vilchis, M.G., Estrada-Soto, S., *et al.* Synthesis of 2-{2-[(α / β -naphthalen-1-ylsulfonyl)amino]-1,3-thiazol-4-yl} acetamides with 11 β -hydroxysteroid dehydrogenase inhibition and in combo antidiabetic activities. *Eur. J. Med. Chem.* **74**, 179-186 (2014).

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, 07/10/2024

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