

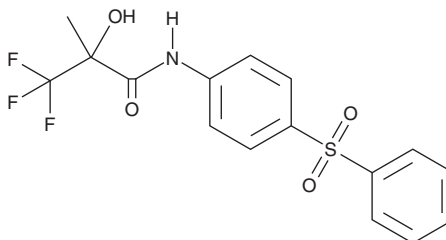
PRODUCT INFORMATION



ZM 226600

Item No. 42440

CAS Registry No.: 147695-92-9
Formal Name: 3,3,3-trifluoro-2-hydroxy-2-methyl-N-[4-(phenylsulfonyl)phenyl]-propanamide
MF: C₁₆H₁₄F₃NO₄S
FW: 373.3
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

ZM 226600 is supplied as a solid. A stock solution may be made by dissolving the ZM 226600 in the solvent of choice, which should be purged with an inert gas. ZM 226600 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

ZM 226600 is an activator of ATP-sensitive potassium channels (K_{ATP}).¹ It induces relaxation of potassium-precontracted isolated guinea pig bladder detrusor muscle strips (EC₅₀ = 0.5 μM). ZM 226600 reduces bladder overactivity in a rat model of ligature-intact partial urethral obstruction.²

References

1. Grant, T.L., Ohnmacht, C.J., and Howe, B.B. Anilide tertiary carbinols: A novel series of K⁺ channel openers. *Trends Pharmacol. Sci.* **15**(11), 402-414 (1994).
2. Pinna, C., Sanvito, P., Bolego, C., *et al.* Effect of the ATP-sensitive potassium channel opener ZM226600 on cystometric parameters in rats with ligature-intact, partial urethral obstruction. *Eur. J. Pharmacol.* **516**(1), 71-77 (2005).

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, 12/09/2024

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