PRODUCT INFORMATION



Acetohexamide

Item No. 23900

CAS Registry No.:	968-81-0	\sim
Formal Name:	4-acetyl-N-[(cyclohexylamino)	
	carbonyl]-benzenesulfonamide	H
MF:	$C_{15}H_{20}N_{2}O_{4}S$	
FW:	324.4	S S
Purity:	≥95%	
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥4 years	0
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

Laboratory Procedures

Acetohexamide is supplied as a solid. A stock solution may be made by dissolving the acetohexamide in the solvent of choice, which should be purged with an inert gas. Acetohexamide is slightly soluble in DMSO.

Description

Acetohexamide is a first generation sulfonylurea that inhibits sulfonylurea receptor 1 (SUR1) linked to the inwardly rectifying potassium channel (K_{IR}6.2) with K_i values of 22.9 and 14.2 μ M in HEK293 cells transfected with the human receptor and in rat brain, respectively.¹ It is metabolized to the hypoglycemic compound L-hydroxyhexamide in vivo.^{2,3} Formulations containing acetohexamide have previously been used in the treatment of type 2 diabetes.

References

- 1. Gopalakrishnan, M., Molinari, E.J., Char-Change, S., et al. Pharmacology of human sulphonylurea receptor SUR1 and inward rectifier K⁺ channel Kir6.2 combination expressed in HEK-293 cells. Br. J. Pharmacol. 129(7), 1323-1332 (2000).
- 2. McMahon, R.E., Marshall, F.J., and Culp, H.W. The nature of the metabolites of acetohexamide in the rat and in the human. J. Pharmacol. Exp. Ther. 149(2), 272-279 (1965).
- 3. Imamura, Y., Sanai, K., Seri, K., et al. Hypoglycemic effect of S(-)-hydroxyhexamide, a major metabolite of acetohexamide, and its enantiomer R(+)-hydroxyhexamide. Life Sci. 69(16), 1947-1955 (2001).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

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

Suyer 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, 11/14/2022

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