

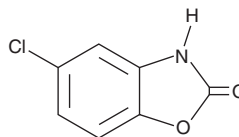
PRODUCT INFORMATION



Chlorzoxazone

Item No. 25826

CAS Registry No.: 95-25-0
Formal Name: 5-chloro-2(3H)-benzoxazolone
Synonym: NSC 26189
MF: C₇H₄ClNO₂
FW: 169.6
Purity: ≥98%
UV/Vis.: λ_{max}: 284 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Chlorzoxazone (Item No. 25826) is an analytical reference standard categorized as a skeletal muscle relaxant.^{1,2} This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 18869).

References

1. Pedarzani, P. and Stocker, M. Molecular and cellular basis of small- and intermediate-conductance, calcium-activated potassium channel function in the brain. *Cell. Mol. Life Sci.* **65(20)**, 3196-3217 (2008).
2. Gao, Z., Todorov, B., Barrett, C.F., *et al.* Cerebellar ataxia by enhanced Ca_v2.1 currents is alleviated by Ca²⁺-dependent K⁺-channel activators in *Cacna1a*^{S218L} mutant mice. *J. Neurosci.* **32(44)**, 15533-15546 (2012).

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, 07/01/2023

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