

PRODUCT INFORMATION



Avizafone (hydrochloride)

Item No. 26527

CAS Registry No.: 60067-16-5

Formal Name: L-lysyl-N-(2-benzoyl-4-chlorophenyl)-N-methyl-glycinamide, dihydrochloride

Synonym: Pro Diazepam

MF: $C_{22}H_{27}ClN_4O_3 \cdot 2HCl$

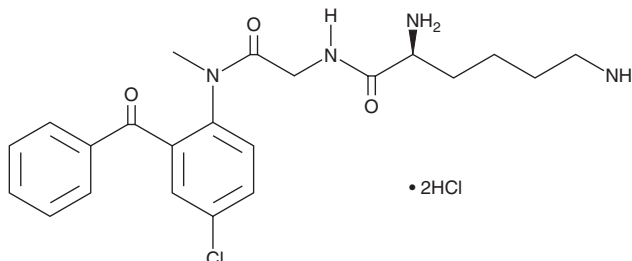
FW: 503.9

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Description

Avizafone (hydrochloride) (Item No. 26527) is an analytical reference standard categorized as a prodrug form of diazepam (Item Nos. ISO60177 | 15887).^{1,2} Formulations containing avizafone have been used as antidotes for nerve agent exposure in a military setting. This product is intended for research and forensic applications.

References

1. Hassall, C.H., Wolmes, S.W., Johnson, W.H., *et al.* Peptido-aminobenzophenones-novel latentiated benzo-1,4-diazepines. *Experientia* **33**(11), 1492-1493 (1977).
2. Price, M.E., Whitmore, C.L., Tattersall, J.E.H., *et al.* Efficacy of the antinicotinic compound MB327 against soman poisoning - Importance of experimental end point. *Toxicol. Lett.* **293**, 167-171 (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, 09/03/2026

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