

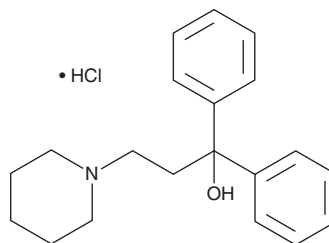
# PRODUCT INFORMATION



## Pridinol (hydrochloride)

Item No. 36056

CAS Registry No.: 968-58-1  
Formal Name:  $\alpha,\alpha$ -diphenyl-1-piperidinepropanol,  
monohydrochloride  
MF:  $C_{20}H_{25}NO \cdot HCl$   
FW: 331.9  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Description

Pridinol (hydrochloride) (Item No. 36056) is an analytical reference standard categorized as a muscle relaxant.<sup>1</sup> Formulations containing pridinol have been used in the treatment of muscle spasms and muscle pain. This product is intended for research and forensic applications.

### Reference

1. Kasuya, Y. Chemicopharmacological studies on antispasmodic action. XII. Structure-activity relationship of aralkylamines. *Chem. Pharm. Bull. (Tokyo)* **6**(2), 147-154 (1958).

#### 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, 07/19/2024

#### 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