

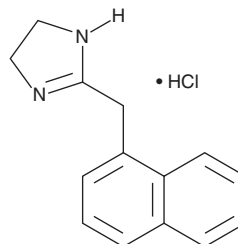
# PRODUCT INFORMATION



## Naphazoline (hydrochloride)

Item No. 23774

**CAS Registry No.:** 550-99-2  
**Formal Name:** 4,5-dihydro-2-(1-naphthalenylmethyl)-1H-imidazole, monohydrochloride  
**Synonym:** 2-(1-Naphthylmethyl)-2-imidazoline  
**MF:**  $C_{14}H_{14}N_2 \cdot HCl$   
**FW:** 246.7  
**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.

### Laboratory Procedures

Naphazoline (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the naphazoline (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Naphazoline (hydrochloride) is slightly soluble in DMSO and methanol.

### Description

Naphazoline is a vasoconstrictor.<sup>1</sup> *In vivo*, naphazoline (0.1 mg/kg, i.v.) decreases specific airway resistance and reverses intranasal blockage induced by U-46619 (Item No. 16450) or leukotriene D<sub>4</sub> (LTD<sub>4</sub>; Item No. 20310) in guinea pigs. Formulations containing naphazoline have been used to treat allergic conjunctivitis and as nasal decongestants.

### Reference

1. Tanaka, Y., Mizutani, N., Fujii, M., *et al.* Different mechanisms between thromboxane A<sub>2</sub>- and leukotriene D<sub>4</sub>-induced nasal blockage in guinea pigs. *Prostaglandins Other Lipid Mediat.* **80(3-4)**, 144-154 (2006).

#### 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, 11/15/2022

#### 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](http://WWW.CAYMANCHEM.COM)