

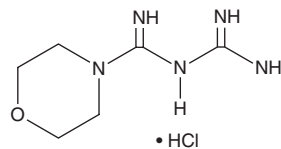
# PRODUCT INFORMATION



## Moroxydine (hydrochloride)

Item No. 21647

**CAS Registry No.:** 3160-91-6  
**Formal Name:** N-(aminoiminomethyl)-4-morpholinecarboximidamide, monohydrochloride  
**Synonym:** ABOB  
**MF:** C<sub>6</sub>H<sub>13</sub>N<sub>5</sub>O • HCl  
**FW:** 207.7  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 241 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Moroxydine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the moroxydine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Moroxydine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of moroxydine (hydrochloride) in these solvents is approximately 0.33, 20, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of moroxydine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of moroxydine (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Moroxydine is a heterocyclic biguanidine compound with antiviral activity. In grass carp reovirus-infected CIK cells, virus-related Bcl-2- and caspase-mediated apoptosis is reduced by 50% following treatment with 6.3 µg/ml of moroxydine for 96 hours.<sup>1</sup> Moroxydine targets DNA and RNA viruses including measles, herpes simplex, varicella zoster, and hepatitis C, and formulations containing moroxydine are used to treat the symptoms of influenza virus strain A infection.<sup>1-3</sup> Moroxydine is considered effective and generally safe in veterinary applications. In gibel carp housed at 25°C, the absorption and elimination half-lives are 2.02 and 4.22 hours, respectively, following oral administration of a 10 mg/kg dose of moroxydine.<sup>4</sup>

### References

1. Yu, X.-B., Chen, X.-H., Ling, F., *et al.* Moroxydine hydrochloride inhibits grass carp reovirus replication and suppresses apoptosis in *Ctenopharyngodon idella* kidney cells. *Antiviral Res.* **131**, 156-165 (2016).
2. Hopkins, E.J., Pye, A.M., Solomon, M., *et al.* Controlled trial of "virugon" in treatment of measles. *Br. Med. J.* **2(5262)**, 1263-1264 (1961).
3. Brown, A.G., Renton, G.H., and Griffiths, P. Comparative study of "Virugon" and penicillin in measles. *Br. Med. J.* **1(5275)**, 374-375 (1962).
4. Liu, W., Xu, J., Zhou, Y., *et al.* Pharmacokinetics and tissue residues of moroxydine hydrochloride in gibel carp, *Carassius gibelio* after oral administration. *J. Vet. Pharmacol. Ther.* **39(4)**, 398-404 (2016).

#### 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, 10/17/2022

#### 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