

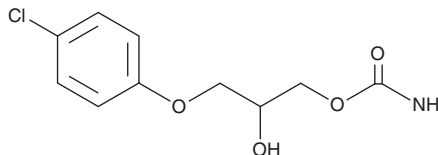
# PRODUCT INFORMATION



## Chlorphenesin Carbamate

Item No. 39252

**CAS Registry No.:** 886-74-8  
**Formal Name:** 3-(4-chlorophenoxy)-1,2-propanediol, 1-carbamate  
**Synonyms:** Maolate, NSC 82943, OC-201  
**MF:** C<sub>10</sub>H<sub>12</sub>ClNO<sub>4</sub>  
**FW:** 245.7  
**Purity:** ≥98%  
**Supplied as:** A 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

Chlorphenesin carbamate is supplied as a solid. A stock solution may be made by dissolving the chlorphenesin carbamate in the solvent of choice, which should be purged with an inert gas. Chlorphenesin carbamate is soluble in methanol.

### Description

Chlorphenesin carbamate is a muscle relaxant.<sup>1</sup> It decreases the dorsal root potential induced by a current-stimulated ventral root, as well as L-glutamate-induced motor neuron discharges, in isolated frog spinal cords when used at a concentration of 1 mM. *In vivo*, chlorphenesin carbamate (50 mg/kg) reduces the amplitude of polysynaptic responses in ipsilateral ventral root L5 induced by electrical stimulation of dorsal root L5 in spinalized rats.<sup>2</sup> It decreases monosynaptic signal transfer from primary afferent fibers to motor neurons in the same animals at the same dose. Formulations containing chlorphenesin carbamate have been used in the treatment of post-surgical pain, acute inflammation, and as a muscle relaxant in veterinary applications.

### References

1. Aihara, H., Kurachi, M., Nakane, S., *et al.* The action of chlorphenesin carbamate on the frog spinal cord. *Jpn. J. Pharmacol.* **30(1)**, 29-36 (1980).
2. Kurachi, M. and Aihara, H. Effect of a muscle relaxant, chlorphenesin carbamate, on the spinal neurons of rats. *Jpn. J. Pharmacol.* **36(1)**, 7-13 (1984).

#### 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, 10/26/2023

#### 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