

PRODUCT INFORMATION



Lercanidipine

Item No. 46189

CAS Registry No.: 100427-26-7

Formal Name: 1,4-dihydro-2,6-dimethyl-4-(3-nitrophenyl)-3,5-pyridinedicarboxylic acid, 3-[2-[(3,3-diphenylpropyl)methylamino]-1,1-dimethylethyl] 5-methyl ester

Synonym: Masnidipine

MF: $C_{36}H_{41}N_3O_6$

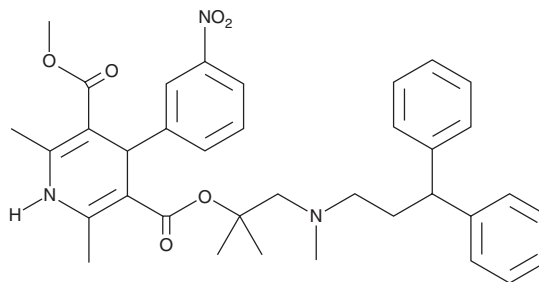
FW: 611.7

Purity: $\geq 98\%$

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

Lercanidipine is supplied as a solid. A stock solution may be made by dissolving the lercanidipine in the solvent of choice, which should be purged with an inert gas. Lercanidipine is soluble (≥ 10 mg/ml) in ethanol and slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Lercanidipine is a dihydropyridine L-type calcium channel blocker.¹ It binds to rat brain and heart homogenates (K_i s = 0.24-0.3 and 0.22 nM, respectively, in radioligand binding assays) and inhibits potassium-induced contraction of isolated rat aorta (IC_{50} = 1.3 nM). Lercanidipine reduces diastolic blood pressure (DBP) in normotensive and spontaneously hypertensive rats with ED_{25} values of 16.3 and 15.5 $\mu\text{g/kg}$, respectively. Formulations containing lercanidipine have been used in the treatment of hypertension.

Reference

1. Testa, R., Leonardi, A., Tajana, A., et al. Lercanidipine (Rec 15/2375): A novel 1,4-dihydropyridine calcium antagonist for hypertension. *Cardiovasc. Drug Rev.* **15**(3), 187-219 (1997).

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, 08/21/2026

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