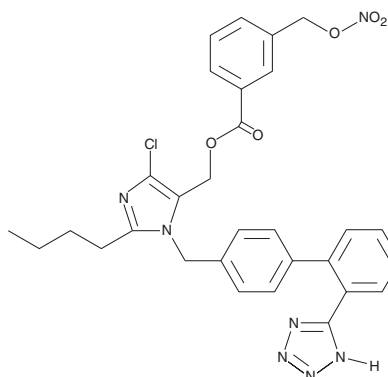


PRODUCT INFORMATION

NO-Losartan A

Item No. 10006456

CAS Registry No.: 791122-48-0
Formal Name: [2-butyl-4-chloro-1-[[2'-(1H-tetrazol-5-yl)[1,1'-biphenyl]-4-yl]methyl]-1H-imidazol-5-yl] methyl ester
MF: C₃₀H₂₈ClN₇O₅
FW: 602.0
Purity: ≥97%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

NO-losartan A is supplied as a crystalline solid. A stock solution may be made by dissolving the NO-losartan A in the solvent of choice, which should be purged with an inert gas. NO-losartan A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NO-losartan A in these solvents is approximately 30 mg/ml.

NO-losartan A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NO-losartan A should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. NO-losartan A has a solubility of approximately 500 µg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Losartan is a mammalian AT₁ receptor antagonist with a K_i value of 5-20 nM.¹ In humans, losartan effectively controls hypertension while protecting renal function.² Nitric oxide (NO) causes vasodilation and also inhibits platelet and neutrophil aggregation in the endothelium.^{3,4} NO-losartan A possesses similar anti-hypertensive effects to losartan, with the addition of the vasodilating effects of NO release.⁵

References

1. Ji, H., Leung, M., Zhang, Y., *et al.* Differential structural requirements for specific binding of nonpeptide and peptide antagonists to the AT₁ angiotensin receptor. Identification of amino acid residues that determine binding of the antihypertensive drug losartan. *J. Biol. Chem.* **269**(24), 16533-16536 (1994).
2. Caruso, D., D'Avino, M., Acampora, C., *et al.* Effects of losartan and chlorthalidone on blood pressure and renal vascular resistance index in non-diabetic patients with essential hypertension and normal renal functions. *J. Cardiovasc. Pharmacol.* **44**(5), 520-524 (2004).
3. Moncada, S., and Higgs, A. The L-arginine-nitric oxide pathway. *N. Engl. J. Med.* **329**, 2002-2012 (1993).
4. Loscalzo, J. and Welch, G. Nitric oxide and its role in the cardiovascular system. *Prog. Cardiovasc. Dis.* **38**, 87-103 (1995).
5. Breschi, M.C., Calderone, V., Digiacomo, M., *et al.* NO-sartans: A new class of pharmacodynamic hybrids as cardiovascular drugs. *J. Med. Chem.* **47**, 5597-5600 (2004).

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, 02/13/2025

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