

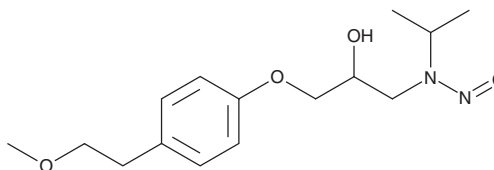
PRODUCT INFORMATION



N-Nitroso Metoprolol

Item No. 34738

CAS Registry No.: 138768-62-4
Formal Name: 1-[4-(2-methoxyethyl)phenoxy]-3-[(1-methylethyl)nitrosoamino]-2-propanol
MF: C₁₅H₂₄N₂O₄
FW: 296.4
Purity: ≥90% (mixture of isomers)
UV/Vis.: λ_{max}: 226 nm
Supplied as: A 10 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

N-Nitroso metoprolol is a genotoxic derivative of metoprolol (Item No. 15429).^{1,2} It induces DNA fragmentation in rat hepatocytes when used at concentrations ranging from 0.1 to 1 mM.¹ *In vivo*, N-nitroso metoprolol (1,000 mg/kg) induces micronucleation of rat hepatocytes but not erythrocytes from the bone marrow or spleen.

References

1. Robbiano, L., Martelli, A., Allavena, A., *et al.* Formation of the N-nitroso derivatives of six β-adrenergic-blocking agents and their genotoxic effects in rat and human hepatocytes. *Cancer Res.* **51(9)**, 2273-2279 (1991).
2. Martelli, A., Allavena, A., Sottofattori, E., *et al.* Low clastogenic activity in vivo of the N-nitroso derivatives of 5 β-adrenergic-blocking drugs proved to be potent genotoxins in vitro. *Toxicol. Lett.* **73(3)**, 185-191 (1994).

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, 03/25/2026

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