

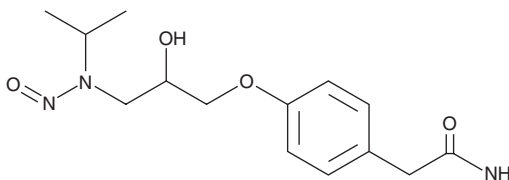
PRODUCT INFORMATION



N-Nitroso Atenolol

Item No. 34735

CAS Registry No.: 134720-04-0
Formal Name: 4-[2-hydroxy-3-[(1-methylethyl)nitrosoamino]propoxy]-benzeneacetamide
Synonym: NO-Atenolol
MF: $C_{14}H_{21}N_3O_4$
FW: 295.3
Purity: $\geq 90\%$
UV/Vis.: λ_{\max} : 229 nm
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

N-Nitroso atenolol is supplied as a solid. A stock solution may be made by dissolving the N-nitroso atenolol in the solvent of choice, which should be purged with an inert gas. N-Nitroso atenolol is soluble in the organic solvent ethanol.

Description

N-Nitroso atenolol is a genotoxic derivative of (\pm)-atenolol (Item No. 17250).¹ It induces DNA fragmentation in rat hepatocytes when used at concentrations ranging from 0.1 to 1 mM.

Reference

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).

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, 12/08/2022

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