

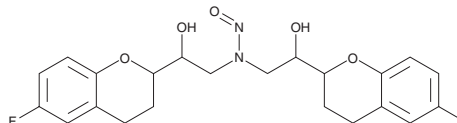
# PRODUCT INFORMATION



## N-Nitroso Nebivolol

Item No. 34739

CAS Registry No.: 1391051-68-5  
Formal Name: N,N-bis(2-(6-fluorochroman-2-yl)-2-hydroxyethyl)nitrous amide  
MF:  $C_{22}H_{24}F_2N_2O_5$   
FW: 434.4  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

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

### Description

N-Nitroso nebivolol is a derivative of the  $\beta_1$ -adrenergic receptor ( $\beta_1$ -AR) antagonist nebivolol (Item No. 23660).

#### 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