

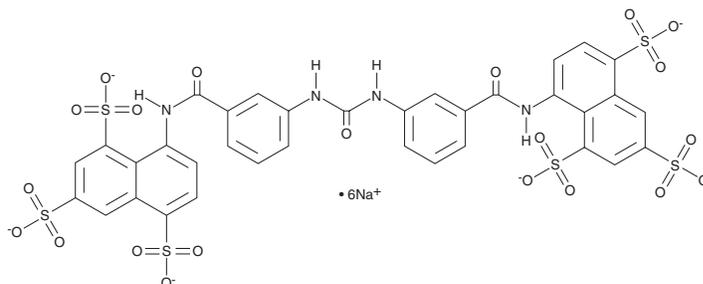
PRODUCT INFORMATION



NF 023

Item No. 17234

CAS Registry No.: 104869-31-0
Formal Name: 8,8'-[carbonylbis(imino-3,1-phenylenecarbonylimino)]bis-1,3,5-naphthalenetrisulfonic acid, hexasodium salt
MF: C₃₅H₂₀N₄O₂₁S₆ • 6Na
FW: 1,162.9
Purity: ≥90%
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

NF 023 is supplied as a solid. A stock solution may be made by dissolving the NF 023 in water. The solubility of NF 023 in water is approximately 100 mM. We do not recommend storing the aqueous solution for more than one day.

Description

NF 023 is a purinergic P2X₁ receptor antagonist (IC₅₀ = 0.21 μM).¹ It is selective for P2X₁ over P2X₂₋₄ receptors (IC₅₀s = >50, 28.9 and >100 μM, respectively). It inhibits α,β-methylene ATP-induced vasoconstriction in isolated rabbit saphenous arteries (pA₂ = 5.69).² NF 023 also inhibits hepatitis C virus (HCV) NS3 helicase (IC₅₀ = 6.2 μM).³

References

1. Soto, F., Lambrecht, G., Nickel, P., *et al.* Antagonistic properties of the suramin analogue NF023 at heterologously expressed P2X receptors. *Neuropharmacology* **38(1)**, 141-149 (1999).
2. Ziyal, R., Ziganshin, A.U., Nickel, P., *et al.* Vasoconstrictor responses via P2X-receptors are selectively antagonized by NF023 in rabbit isolated aorta and saphenous artery. *Br. J. Pharmacol.* **120(5)**, 954-960 (1997).
3. Mukherjee, S., Hanson, A.M., Shadrack, W.R., *et al.* Identification and analysis of hepatitis C virus NS3 helicase inhibitors using nucleic acid binding assays. *Nucleic Acids Res.* **40(17)**, 8607-8621 (2012).

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, 11/15/2022

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