

PRODUCT INFORMATION



NS 3728

Item No. 39988

CAS Registry No.: 265646-85-3

Formal Name: N-[3,5-bis(trifluoromethyl)phenyl]-
N'-[4-bromo-2-(2H-tetrazol-5-yl)
phenyl]-urea

Synonyms: Endovion, SCO 101

MF: C₁₆H₉BrF₆N₆O

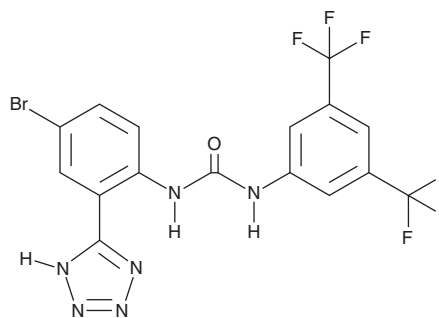
FW: 495.2

Purity: ≥95%

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

NS 3728 is supplied as a solid. A stock solution may be made by dissolving the NS 3728 in the solvent of choice, which should be purged with an inert gas. NS 3728 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

NS 3728 is a chloride channel inhibitor.^{1,2} It inhibits the volume-regulated anion channel (VRAC) in HEK293 cells and chloride conductance in washed isolated human erythrocytes (IC₅₀s = 0.4 and 0.6 μM, respectively).¹ NS 3728 also inhibits calcium-activated chloride currents in Ehrlich-Lettre ascites carcinoma cells (IC₅₀s = 2.11 and 0.7 μM at +95 and -50 mV, respectively).² It reduces the proliferation of Ehrlich-Lettre ascites cells and inhibits IFN-γ-induced nitric oxide (NO) production in BV-2 microglia (IC₅₀s = 0.52 and 3.7 μM, respectively).^{2,3}

References

1. Hélix, D., Strøbaek, B.H.D., and Christophersen, P. Inhibition of the endogenous volume-regulated anion channel (VRAC) in HEK293 cells by acidic di-aryl-ureas. *J. Membr. Biol.* **196**(2), 83-94 (2003).
2. Klausen, T.K., Bergdahl, A., Hougaard, C., *et al.* Cell cycle-dependent activity of the volume- and Ca²⁺-activated anion currents in Ehrlich Lettre ascites cells. *J. Cell. Physiol.* **210**(3), 831-842 (2007).
3. Kjaer, K., Strøbaek, D., Christophersen, P., *et al.* Chloride channel blockers inhibit iNOS expression and NO production in IFNγ-stimulated microglial BV2 cells. *Brain Res.* **1281**, 15-24 (2009).

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, 03/20/2024

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