

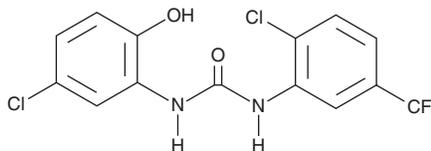
PRODUCT INFORMATION



NS 1738

Item No. 21018

CAS Registry No.: 501684-93-1
Formal Name: N-(5-chloro-2-hydroxyphenyl)-N'-[2-chloro-5-(trifluoromethyl)phenyl]-urea
Synonym: NSC 213859
MF: C₁₄H₉Cl₂F₃N₂O₂
FW: 365.1
Purity: ≥98%
UV/Vis.: λ_{max}: 209, 263, 296 nm
Supplied as: A crystalline 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 1738 is supplied as a crystalline solid. A stock solution may be made by dissolving the NS 1738 in the solvent of choice, which should be purged with an inert gas. NS 1738 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of NS 1738 in ethanol is approximately 10 mg/ml and approximately 25 mg/ml in DMSO and DMF.

NS 1738 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NS 1738 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. NS 1738 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NS 1738 is a positive allosteric modulator of the α7-containing neuronal nicotinic acetylcholine receptors (nAChRs).¹ When applied in the presence of ACh, NS 1738 increases the peak amplitude of the current flowing through α7-containing nAChRs (EC₅₀ = 3.4 μM).¹ NS 1738 displays no substantial activity for α4β2-, α3β3-, and α1-containing receptors. It is modestly brain-penetrant and counteracts (-)-scopolamine-induced deficit in acquisition of a water-maze learning task in rats.¹ NS 1738 also improves performance in the rat social recognition test.¹ NS 1738 is used to selectively evaluate the role of α7-containing nAChRs in signaling pathways.^{2,3}

References

1. Timmermann, D.B., Gronlien, J.H., Kohlhaas, K.L., *et al.* An allosteric modulator of the α7 nicotinic acetylcholine receptor possessing cognition-enhancing properties *in vivo*. *JPET* **323**(1), 294-307 (2007).
2. Cheng, Q. and Yakel, J.L. Presynaptic α7 nicotinic acetylcholine receptors enhance hippocampal mossy fiber glutamatergic transmission via PKA activation. *J. Neurosci.* **34**(1), 124-133 (2014).
3. Freitas, K., Carroll, F.I., and Damaj, M.I. The antinociceptive effects of nicotinic receptors α7-positive allosteric modulators in murine acute and tonic pain models. *J. Pharmacol. Exp. Ther.* **344**(1), 264-275 (2013).

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, 11/14/2022

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