

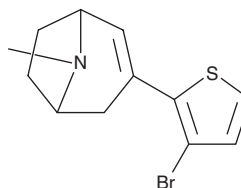
PRODUCT INFORMATION



NS 3861

Item No. 37406

CAS Registry No.: 216853-59-7
Formal Name: 3-(3-bromo-2-thienyl)-8-methyl-8-azabicyclo[3.2.1]oct-2-ene
MF: $C_{12}H_{14}BrNS$
FW: 284.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

NS 3861 is supplied as a solid. A stock solution may be made by dissolving the NS 3861 in the solvent of choice, which should be purged with an inert gas. NS 3861 is soluble in methanol, ethanol, and DMSO.

Description

NS 3861 is a neuronal nicotinic acetylcholine receptor (nAChR) agonist.¹ It induces inward currents in patch-clamp assays using *X. laevis* oocytes expressing recombinant human $\alpha 3\beta 2$ - or $\alpha 3\beta 4$ subunit-containing nAChR receptors ($EC_{50}s = 1.6$ and $1 \mu M$, respectively). *In vivo*, NS 3861 (0.1 mg/kg) increases the number of fecal pellets expelled in mice exposed to long-term, but not short-term, morphine injections.²

References

1. Harpsøe, K., Hald, H., Timmermann, D.B., *et al.* Molecular determinants of subtype-selective efficacies of cytosine and the novel compound NS3861 at heteromeric nicotinic acetylcholine receptors. *J. Biol. Chem.* **288**(4), 2559-2570 (2013).
2. Gade, A.R., Kang, M., Khan, F., *et al.* Enhanced sensitivity of $\alpha 3\beta 4$ nicotinic receptors in enteric neurons after long-term morphine: Implication for opioid-induced constipation. *J. Pharmacol. Exp. Ther.* **357**(3), 520-528 (2016).

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, 11/07/2023

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