

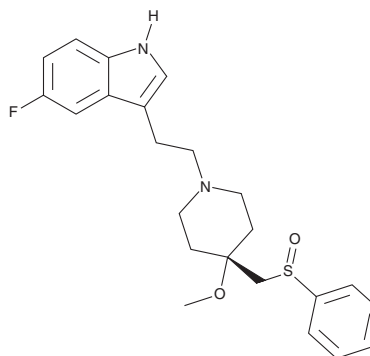
PRODUCT INFORMATION



GR159897

Item No. 32774

CAS Registry No.: 158848-32-9
Formal Name: 5-fluoro-3-[2-[4-methoxy-4-[[[*(R)*-phenylsulfinyl]methyl]-1-piperidinyl]ethyl]-1H-indole
MF: C₂₃H₂₇FN₂O₂S
FW: 414.5
Purity: ≥98%
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

GR159897 is supplied as a solid. A stock solution may be made by dissolving the GR159897 in the solvent of choice, which should be purged with an inert gas. GR159897 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

GR159897 is a nonpeptide antagonist of the neurokinin-2 (NK₂) receptor.¹ It binds to NK₂ receptors (K_i = 0.32 nM) and is selective for NK₂ over NK₁ and NK₃ receptors in radioligand binding assays (K_is = 5,010 and >1,000 nM, respectively). GR159897 inhibits contractions induced by the NK₂ receptor agonist GR64349 in isolated guinea pig trachea (pA₂ = 8.7). *In vivo* GR159897 (0.12 mg/kg) inhibits GR64349-induced bronchoconstriction in anesthetized guinea pigs. GR159897 (0.2-50 µg/kg) increases the amount of time spent in the front of the cage during confrontation with a human observer, a marker of anxiolytic-like activity, in marmosets.²

References

1. Beresford, I.J., Sheldrick, R.L., Ball, D.I., *et al.* GR159897, a potent non-peptide antagonist at tachykinin NK₂ receptors. *Eur. J. Pharmacol.* **272(2-3)**, 241-248 (1995).
2. Walsh, D.M., Stratton, S.C., Harvey, F.J., *et al.* The anxiolytic-like activity of GR159897, a non-peptide NK₂ receptor antagonist, in rodent and primate models of anxiety. *Psychopharmacol. (Berl)*. **121(2)**, 186-191 (1995).

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, 12/13/2022

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