

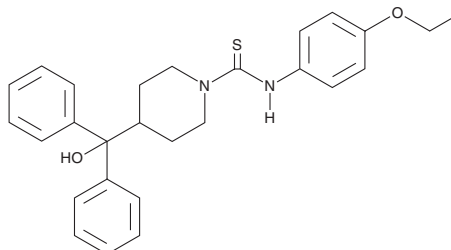
PRODUCT INFORMATION



SF11

Item No. 33355

CAS Registry No.: 443292-81-7
Formal Name: N-(4-ethoxyphenyl)-4-(hydroxydiphenylmethyl)-1-piperidinecarbothioamide
MF: C₂₇H₃₀N₂O₂S
FW: 446.6
Purity: ≥95%
UV/Vis.: λ_{max}: 252 nm
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

SF11 is supplied as a solid. A stock solution may be made by dissolving the SF11 in the solvent of choice, which should be purged with an inert gas. SF11 is soluble in DMSO.

Description

SF11 is an antagonist of the neuropeptide Y (NPY) receptor Y₂ (IC₅₀ = 0.199 μM).¹ It is selective for Y₂ over Y₁ receptors (IC₅₀ = >33 μM), as well as a panel of 40 additional receptors, ion channels, and transporters at 10 μM. SF11 (10 and 30 mg/kg) decreases immobility time in the forced swim test in rats.²

References

1. Brothers, S.P., Saldanha, A., Spicer, T.P., *et al.* Selective and brain penetrant neuropeptide Y Y2 receptor antagonists discovered by whole-cell high-throughput screening. *Mol. Pharmacol.* **77**(1), 46-57 (2010).
2. Domin, H., Piergies, N., Pięta, E., *et al.* Characterization of the brain penetrant neuropeptide Y Y2 receptor antagonist SF-11. *ACS Chem. Neurosci.* **10**(8), 3454-3463 (2019).

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