

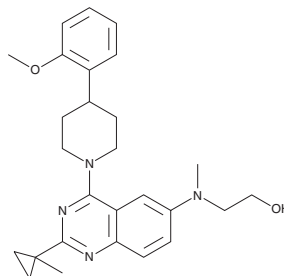
PRODUCT INFORMATION



SBI-810

Item No. 45376

CAS Registry No.: 1849603-79-7
Formal Name: 2-[[4-[4-(2-methoxyphenyl)-1-piperidinyl]-2-(1-methylcyclopropyl)-6-quinazoliny]methylamino]-ethanol
MF: C₂₇H₃₄N₄O₂
FW: 446.6
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

SBI-810 is supplied as a solid. A stock solution may be made by dissolving the SBI-810 in the solvent of choice, which should be purged with an inert gas. SBI-810 is slightly soluble (0.1-1 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

Description

SBI-810 is a β -arrestin-2-biased positive allosteric modulator of neurotensin receptor 1 (NTS₁).^{1,2} It increases neurotensin-1-induced β -arrestin-2 recruitment to human NTS₁ receptors in HEK293T cells (EC₅₀ = 98 nM).² SBI-810 (12 mg/kg) increases the paw withdrawal threshold in mouse models of postoperative, neuropathic, and inflammatory pain.¹ It reduces morphine-induced conditioned place preference (CPP) in mice without inducing CPP itself when administered at a dose of 30 mg/kg.

References

1. Guo, R., Chen, O., Zhou, Y., *et al.* Arrestin-biased allosteric modulator of neurotensin receptor 1 alleviates acute and chronic pain. *Cell* **188(16)**, 4332-4349.e21 (2025).
2. Pinkerton, A.B., Peddibhotla, S., Yamamoto, F., *et al.* Discovery of β -arrestin biased, orally bioavailable, and CNS penetrant neurotensin receptor 1 (NTR1) allosteric modulators. *J. Med. Chem.* **62(17)**, 8357-8363 (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, 04/02/2026

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