

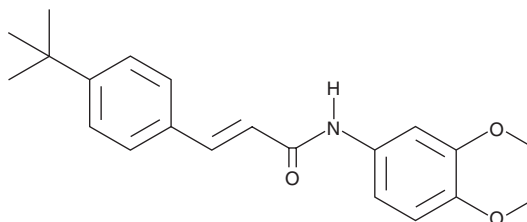
PRODUCT INFORMATION



AMG 9810

Item No. 14715

CAS Registry No.: 545395-94-6
Formal Name: (2E)-N-(2,3-dihydro-1,4-benzodioxin-6-yl)-3-[4-(1,1-dimethylethyl)phenyl]-2-propenamamide
MF: C₂₁H₂₃NO₃
FW: 337.4
Purity: ≥98%
UV/Vis.: λ_{max}: 304 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

AMG 9810 is supplied as a crystalline solid. A stock solution may be made by dissolving the AMG 9810 in the solvent of choice, which should be purged with an inert gas. AMG 9810 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AMG 9810 in ethanol is approximately 10 mg/ml and approximately 16 mg/ml in DMSO and DMF.

Description

AMG 9810 is a competitive antagonist of capsaicin activation of the vanilloid receptor 1 (TRPV1) (IC₅₀s = 24.5 and 85.6 nM for human and rat, respectively).¹ Additionally, it has been shown to compete with other modes of TRPV1 activation, including protons (IC₅₀s = 92.7 and 294 nM for human and rat, respectively), heat (IC₅₀s = 15.8 and 21 nM for human and rat, respectively), and the endogenous ligands: anandamide, N-arachidonyl dopamine (IC₅₀s = 8.5 and 260 nM for human and rat, respectively), and oleoyldopamine.¹ In a rat model of inflammatory pain induced by intraplantar injection of complete Freund's adjuvant, 30-100 mg/kg AMG 9810 reverses thermal and mechanical hyperalgesia.¹

Reference

1. Gavva, N.R., Tamir, R., Qu, Y., *et al.* AMG 9810 [(E)-3-(4-t-butylphenyl)-N-(2,3-dihydrobenzo[b][1,4]dioxin-6-yl)acrylamide], a novel vanilloid receptor 1 (TRPV1) antagonist with antihyperalgesic properties. *J. Pharmacol. Exp. Ther.* **313**(1), 474-484 (2005).

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, 10/31/2022

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