

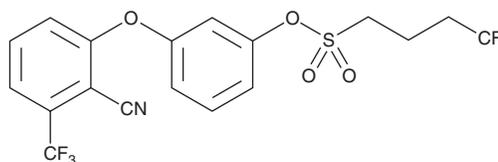
PRODUCT INFORMATION



BAY 59-3074

Item No. 21036

CAS Registry No.: 406205-74-1
Formal Name: 4,4,4-trifluoro-1-butananesulfonic acid, 3-[2-cyano-3-(trifluoromethyl)phenoxy]phenyl ester
MF: C₁₈H₁₃F₆NO₄S
FW: 453.4
Purity: ≥98%
UV/Vis.: λ_{max}: 294 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

BAY 59-3074 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY 59-3074 in the solvent of choice, which should be purged with an inert gas. BAY 59-3074 is soluble in organic solvents such as ethanol and DMSO. The solubility of BAY 59-3074 in these solvents is approximately 100 mM.

Description

BAY 59-3074 is a partial agonist of the cannabinoid (CB) receptors CB₁ and CB₂ (K_is = 55.4, 48.3, and 45.5 nM for rat CB₁, human CB₁, and human CB₂, respectively).¹ BAY 59-3074 is orally active and exhibits anti-hyperalgesic and anti-allodynic properties in rat models of chronic neuropathic and inflammatory pain.¹⁻³

References

1. De Vry, J. and Jentzsch, K.R. Discriminative stimulus effects of the structurally novel cannabinoid CB₁/CB₂ receptor partial agonist BAY 59-3074 in the rat. *Eur. J. Pharmacol.* **505(1-3)**, 127-133 (2004).
2. Teng, H., Thakur, G.A., and Makriyannis, A. Conformationally constrained analogs of BAY 59-3074 as novel cannabinoid receptor ligands. *Bioorg. Med. Chem. Lett.* **21(19)**, 5999-6002 (2011).
3. De Vry, J., Denzer, D., Reissmueller, E., et al. 3-[2-Cyano-3-(trifluoromethyl)phenoxy]phenyl-4,4,4-trifluoro-1-butananesulfonate (BAY 59-3074): A novel cannabinoid CB₁/CB₂ receptor partial agonist with antihyperalgesic and antiallodynic effects. *J. Pharmacol. Exp. Ther.* **310(2)**, 620-632 (2004).

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, 10/05/2022

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