

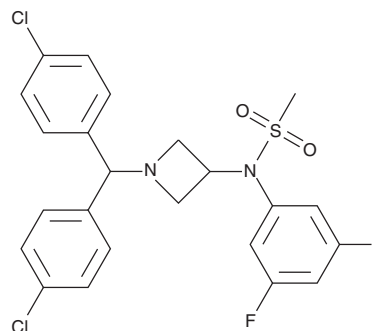
PRODUCT INFORMATION



AVE-1625

Item No. 10009021

CAS Registry No.: 358970-97-5
Formal Name: N-[1-[bis(4-chlorophenyl)methyl]-3-azetidiny]-N-(3,5-difluorophenyl)methanesulfonamide
Synonym: Drinabant
MF: C₂₃H₂₀Cl₂F₂N₂O₂S
FW: 497.4
Purity: ≥98%
UV/Vis.: λ_{max}: 231 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

AVE-1625 is supplied as a crystalline solid. A stock solution may be made by dissolving the AVE-1625 in the solvent of choice, which should be purged with an inert gas. AVE-1625 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AVE-1625 in ethanol is approximately 0.15 mg/ml and approximately 15 mg/ml in DMSO and DMF.

AVE-1625 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AVE-1625 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AVE-1625 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The central cannabinoid (CB₁) receptor is a G protein-coupled receptor that is widely distributed in the central nervous system and several peripheral tissues and binds the active component of *cannabis*, Δ⁹-tetrahydrocannabinol.¹ Signaling through the CB₁ receptor is implicated in attentional and working memory deficits as well as obesity.²⁻⁴ AVE-1625 is a highly potent, selective antagonist for the CB₁ receptor with K_i values of 0.16-0.44 nM.⁴ At 1-3 mg/kg, AVE-1625 significantly improves the performance of rodents in working memory tasks.⁴ At 30 mg/kg, AVE-1625 reduces caloric intake by more than 50% of controls and significantly increases lipolysis from fat tissues and reduces hepatic glycogen levels in rodents.⁵

References

1. Howlett, A.C., Song, C., Berglund, B.A., *et al.* Characterization of CB₁ cannabinoid receptors using receptor peptide fragments and site-directed antibodies. *Mol. Pharmacol.* **53**(3), 504-510 (1998).
2. Williams, J.L., Kashfi, K., Ouyang, N., *et al.* NO-donating aspirin inhibits intestinal carcinogenesis in min (APC^{Min/+}) mice. *Biochem. Biophys. Res. Commun.* **313**(3), 784-788 (2004).
3. Mackie, K. Cannabinoid receptors as therapeutic targets. *Annu. Rev. Pharmacol. Toxicol.* **46**, 101-122 (2006).
4. Borowsky, B., Stevens, R., Mark, B., *et al.* AVE1625, a cannabinoid CB₁ antagonist, as a co-treatment for Schizophrenia: Improvement in cognitive function and reduction of antipsychotic-side effects in animal models. *Neuropsychopharmacology* **30**, S116-S117 (2005).
5. Herling, A.W., Gossel, M., Haschke, G., *et al.* The CB₁ receptor antagonist AVE1625 affects primarily metabolic parameters independently of reduced food intake in wistar rats. *Am. J. Physiol. Endocrinol. Metab.* **293**(3), E826-E832 (2007).

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, 01/24/2024

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