

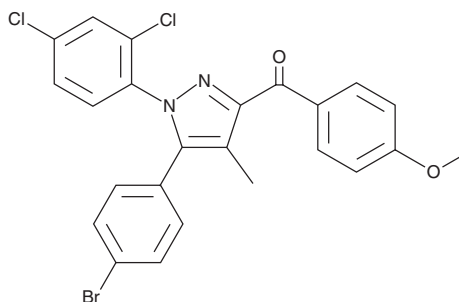
PRODUCT INFORMATION



ABD459

Item No. 20003

CAS Registry No.: 1047670-51-8
Formal Name: [5-(4-bromophenyl)-1-(2,4-dichlorophenyl)-4-methyl-1H-pyrazol-3-yl](4-methoxyphenyl)methanone
MF: C₂₄H₁₇BrCl₂N₂O₂
FW: 516.2
Purity: ≥95%
UV/Vis.: λ_{max}: 249, 295 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

ABD459 is supplied as a crystalline solid. A stock solution may be made by dissolving the ABD459 in the solvent of choice. ABD459 is soluble in the organic solvent dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ABD459 in DMF is approximately 3 mg/ml.

ABD459 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ABD459 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ABD459 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ABD459 is a neutral antagonist of the central cannabinoid 1 (CB₁) receptor (K_i = 8.6 nM).¹ It inhibits food consumption in nonfasted mice without affecting motor activity.¹ ABD459 reduces active food seeking for 5-6 hours after treatment, with no rebound after washout. ABD459 also diminishes rapid eye movement (REM) sleep, with no alterations of wakefulness or non-REM sleep.¹

Reference

1. Goonawardena, A. V., Plano, A. Robinson, L., *et al.* Modulation of food consumption and sleep-wake in mice by the neutral CB1 antagonist ABD459. *Behav. Pharmacol.* **26**(3), 289-303 (2015).

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, 02/15/2024

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