

PRODUCT INFORMATION



JJH260

Item No. 19384

CAS Registry No.: 1831135-30-8
Formal Name: 4-[2-(4-chlorophenyl)ethyl]-1-piperidinecarboxylic acid, 7-[4-(dimethylamino)benzoyl]hexahydro-1,3-dioxoimidazo[1,5-a]pyrazin-2-yl ester

MF: C₂₉H₃₄ClN₅O₅

FW: 568.1

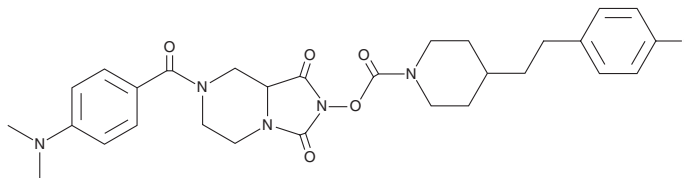
Purity: ≥90%

UV/Vis.: λ_{max}: 220, 301 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

JJH260 is supplied as a crystalline solid. A stock solution may be made by dissolving the JJH260 in the solvent of choice. JJH260 is soluble in the organic solvent dimethyl formamide (DMF), which should be purged with an inert gas, at a concentration of approximately 1 mg/ml.

JJH260 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JJH260 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JJH260 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

JJH260 is an N-hydroxy hydantoin carbamate that inhibits androgen-induced gene 1 (AIG1), an enzyme that hydrolyzes fatty acid esters of hydroxy fatty acids (FAHFAs). It blocks the hydrolysis of 9-PAHSA (Item No. 17037) with an IC₅₀ value of 0.57 μM.¹ JJH260 also inhibits the novel FAHFA hydrolase androgen-dependent TFPI-regulating protein (ADTRP; IC₅₀ = 8.5 μM), as well as the serine hydrolase ABHD6 and the lysophospholipases LYPLA1 and LYPLA2.¹ JJH260 inhibits the FAHFA hydrolase activity of LNCaP and T cell lysates and intact cells.¹

Reference

1. Parsons, W.H., Kolar, M.J., Kamat, S.S., *et al.* AIG1 and ADTRP are atypical integral membrane hydrolases that degrade bioactive FAHFAs. *Nat. Chem. Biol.* **12**(5), 367-372 (2016).

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