

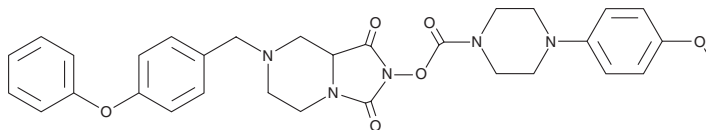
PRODUCT INFORMATION



ABC34

Item No. 19385

CAS Registry No.: 1831135-56-8
Formal Name: 4-(4-methoxyphenyl)-1-piperazinecarboxylic acid, hexahydro-1,3-dioxo-7-[[4-phenoxyphenyl)methyl]imidazo[1,5-a]pyrazin-2-yl ester
MF: C₃₁H₃₃N₅O₆
FW: 571.6
Purity: ≥98%
UV/Vis.: λ_{max}: 235 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

ABC34 is supplied as a crystalline solid. A stock solution may be made by dissolving the ABC34 in the solvent of choice. ABC34 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ABC34 in these solvents is approximately 10 and 20 mg/ml, respectively.

ABC34 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ABC34 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ABC34 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ABC34 is an inactive control probe for JJH260 (Item No. 19384), the inhibitor of androgen-induced gene 1 (AIG1), an enzyme that hydrolyzes fatty acid esters of hydroxy fatty acids (FAHFAs).¹ ABC34 demonstrates an IC₅₀ value greater than 25 μM for the inhibition of 9-PAHSA (Item No. 17037) but is more potent in blocking the serine hydrolase ABHD6 and the palmitoyl-protein hydrolase PPT1, which are both off-targets of JJH260.¹

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.* **3.28**, (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