

PRODUCT INFORMATION



JP104

Item No. 10008661

CAS Registry No.: 887264-45-1

Formal Name: N-10-undecyn-1-yl-carbamic acid, 3'-(aminocarbonyl)[1,1'-biphenyl]-3-yl ester

Synonym: Click Tag™ JP104

MF: C₂₅H₃₀N₂O₃

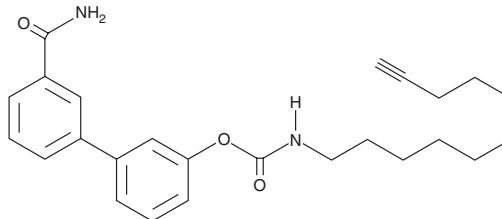
FW: 406.5

Purity: ≥98%

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

JP104 is supplied as a crystalline solid. A stock solution may be made by dissolving the JP104 in an organic solvent purged with an inert gas. JP104 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JP104 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

JP104 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JP104 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. JP104 has a solubility of approximately 0.25 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 enzyme, fatty acid amide hydrolase (FAAH), is widely expressed in brain and other tissues, and is capable of hydrolyzing anandamide (AEA) and other simple esters and amides with long unsaturated acyl chains.¹ JP104 is an irreversible FAAH inhibitor of the carbamate class with an IC₅₀ of 7.3 nM for the human recombinant enzyme when tested using radiolabeled oleamide as the substrate.² The alkyl derivative on JP104 reacts with azide-modified reporter tags, such as azido-rhodamine or azido-biotin, for visualization of JP104 bound to FAAH *in vivo*.

References

1. Zhang, Y.H., Zhang, G.Y., Mollat, P., *et al.* Purification, characterization, and cellular localization of the 100-kDa human placental GTPase-activating protein. *J. Biol. Chem.* **268**, 18875-18881 (1993).
2. Alexander, J.P. and Cravatt, B.F. Mechanism of carbamate inactivation of FAAH: Implications for the design of covalent inhibitors and *in vivo* functional probes for enzymes. *Chemistry & Biology* **12**, 1179-1187 (2005).

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, 05/01/2024

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