

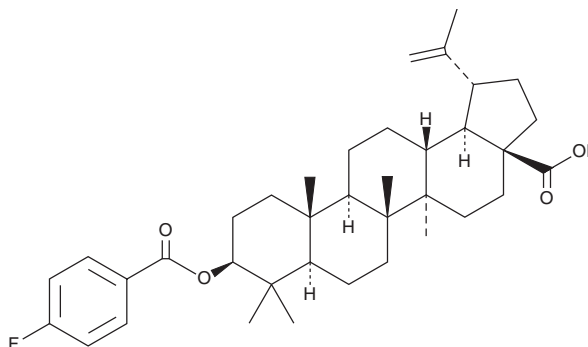
PRODUCT INFORMATION



ZM 522

Item No. 37904

CAS Registry No.: 1454575-38-2
Formal Name: 3β-[(4-fluorobenzoyl)oxy]-lup-20(29)-en-28-oic acid
MF: C₃₇H₅₁FO₄
FW: 578.8
Purity: ≥98%
UV/Vis.: λ_{max}: 241 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

ZM 522 is supplied as a solid. A stock solution may be made by dissolving the ZM 522 in the solvent of choice, which should be purged with an inert gas. ZM 522 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ZM 522 in ethanol and DMSO is approximately 2 mg/ml and approximately 10 mg/ml in DMF.

Description

ZM 522 is an inhibitor of ecto-5'-nucleotidase, also known as CD73 (IC₅₀s = 0.56 and 3.06 μM for the human and mouse enzymes, respectively), and a derivative of the triterpenoid betulinic acid (Item No. 11686).¹ It increases the levels of IFN-γ secreted by isolated mouse spleen cells when used at a concentration of 10 μM in the presence of AMP.

Reference

1. Zhang, Y., Yang, K., Ye, S., *et al.* Application of a fluorine strategy in the lead optimization of betulinic acid to the discovery of potent CD73 inhibitors. *Steroids* **188**, 109112 (2022).

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, 04/26/2023

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