

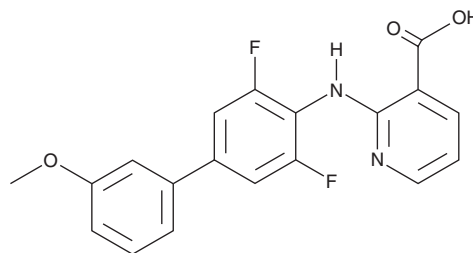
PRODUCT INFORMATION



ASLAN003

Item No. 33516

CAS Registry No.: 1035688-66-4
Formal Name: 2-[(3,5-difluoro-3'-methoxy[1,1'-biphenyl]-4-yl)amino]-3-pyridinecarboxylic acid
MF: C₁₉H₁₄F₂N₂O₃
FW: 356.3
Purity: ≥98%
UV/Vis.: λ_{max}: 288 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

ASLAN003 is supplied as a crystalline solid. A stock solution may be made by dissolving the ASLAN003 in the solvent of choice, which should be purged with an inert gas. ASLAN003 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ASLAN003 in these solvents is approximately 30 mg/ml.

ASLAN003 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ASLAN003 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ASLAN003 has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ASLAN003 is an inhibitor of dihydroorotate dehydrogenase (DHODH; IC₅₀ = 35 nM).¹ It inhibits the proliferation of MOLM-14 and KG-1 acute myeloid leukemia (AML) cells (IC₅₀s = 0.582 and 0.382 μM, respectively), as well as induces apoptosis in the same cells when used at concentrations of 0.5 and 1 μM. ASLAN003 (1, 2, and 4 μM) induces differentiation in bone marrow cells isolated from patients with AML or myelodysplastic syndrome (MDS). It improves survival in a patient-derived xenograft (PDX) mouse model of AML when administered at a dose of 50 mg/kg.

Reference

1. Zhou, J., Quah, J.Y., Ng, Y., *et al.* ASLAN003, a potent dihydroorotate dehydrogenase inhibitor for differentiation of acute myeloid leukemia. *Haematologica* **105(9)**, 2286-2297 (2020).

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, 10/31/2022

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