

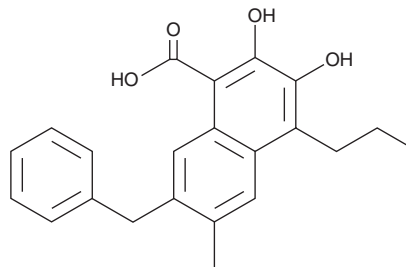
PRODUCT INFORMATION



FX11

Item No. 38518

CAS Registry No.: 213971-34-7
Formal Name: 2,3-dihydroxy-6-methyl-7-(phenylmethyl)-4-propyl-1-naphthalenecarboxylic acid
MF: C₂₂H₂₂O₄
FW: 350.4
Purity: ≥95%
UV/Vis.: λ_{max}: 233 nm
Supplied as: A 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

FX11 is supplied as a solid. A stock solution may be made by dissolving the FX11 in the solvent of choice, which should be purged with an inert gas. FX11 is soluble in DMSO.

Description

FX11 is an inhibitor of lactate dehydrogenase-A (LDH-A; $K_i = 8 \mu\text{M}$).¹ It is selective for LDH-A over GAPDH ($K_i = >300 \mu\text{M}$). FX11 increases oxygen consumption and the production of reactive oxygen species (ROS), as well as induces cell death in P493-6 lymphoblastoid cells. *In vivo*, FX11 (42 $\mu\text{g}/\text{animal}$, i.p.) reduces tumor volume in a P493-6 mouse xenograft model. It also reduces tumor weight in a BxPC-3 mouse xenograft model when administered at a dose of 2 mg/kg.²

References

1. Le, A., Cooper, C.R., Gouw, A.M., *et al.* Inhibition of lactate dehydrogenase A induces oxidative stress and inhibits tumor progression. *Proc. Nat. Acad. Sci. USA* **107**(5), 2037-2047 (2009).
2. Mohammed, G.H., Vassileva, V., Acedo, P., *et al.* Targeting pyruvate kinase M2 and lactate dehydrogenase A is an effective combination strategy for the treatment of pancreatic cancer. *Cancers (Basel)* **11**(9), 1372 (2019).

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, 07/26/2023

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