

YM-53601

YM-53601 inhibits squalene synthase activities

YM-53601, a squalene synthase inhibitor, reduces plasma cholesterol and triglyceride levels in vivo. YM-53601 inhibits squalene synthase derived from human hepatoma cells. Lipid-lowering agent. YM-53601 is also an inhibitor of farnesyl-diphosphate farnesyltransferase 1 (FDFT1) enzyme activity and abrogates HCV propagation.

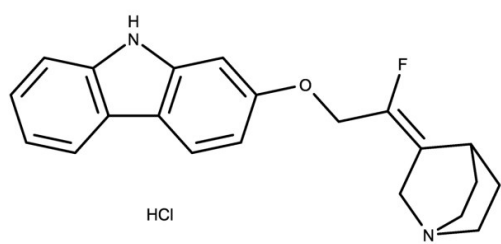
Ordering Information

[Order Online »](#)

ENZ-CHM463-0001	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 9H-Carbazole, 2-[(2E)-2-(1-azabicyclo[2.2.2]oct-3-ylidene)-2-fluoroethoxy]-, hydrochloride (1:1)

Appearance White to light brown solid.

CAS 182959-33-7

Couple Target Squalene synthase

Couple Type Inhibitor

Formula $C_{21}H_{22}ClFN_2O$

Identity Determined by NMR.

MW 372.86

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (100mg/ml).

Last modified: January 30, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)