

Emricasan

Pan-Caspase inhibitor

Emricasan is a potent, selective, and orally bioavailable pan-caspase inhibitor. It is designed to inhibit the activity of caspases, which are enzymes involved in the process of apoptosis (programmed cell death) and inflammation.

Emricasan has been investigated for its potential therapeutic applications in several conditions, including:

- **Liver Diseases:** Chronic liver diseases such as non-alcoholic steatohepatitis (NASH), liver fibrosis, and cirrhosis.
- **Cardiovascular Diseases:** Conditions involving cardiac apoptosis and inflammation.
- **Neurodegenerative Diseases:** Diseases characterized by excessive neuronal apoptosis, such as Alzheimer's disease and Huntington's disease.

Ordering Information

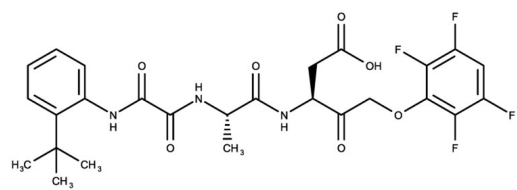
[Order Online »](#)

ENZ-CHM348-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- GMP format available



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20°C for up to 1 month.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(3S)-3-[[[(2S)-2-[[2-(2-tert-butylanilino)-2-oxoacetyl]amino]propanoyl]amino]-4-oxo-5-(2,3,5,6-tetrafluorophenoxy)pentanoic acid, PF 03491390, IDN-6556
Appearance	White solid.
CAS	254750-02-2
Couple Target	Caspase
Couple Type	Inhibitor
Formula	$C_{26}H_{27}F_4N_3O_7$
Identity	Determined by NMR.
MW	569.51
Purity	≥98% HPLC
Solubility	Soluble in DMSO (20 mg/ml).



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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)