

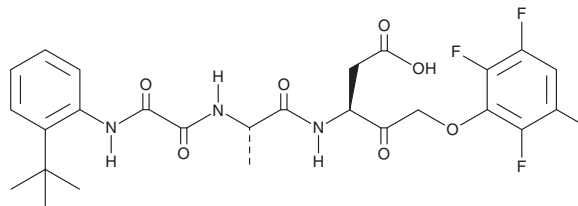
PRODUCT INFORMATION



Emricasan

Item No. 22204

CAS Registry No.: 254750-02-2
Formal Name: N-[2-(1,1-dimethylethyl)phenyl]-2-oxoglycyl-N-[(1S)-1-(carboxymethyl)-2-oxo-3-(2,3,5,6-tetrafluorophenoxy)propyl]-L-alaninamide
Synonyms: IDN-6556, PF-03491390
MF: C₂₆H₂₇F₄N₃O₇
FW: 569.5
Purity: ≥95%
UV/Vis.: λ_{max}: 266 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

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

Description

Emricasan is a pan-caspase inhibitor.¹⁻⁴ *Ex vivo*, emricasan (10 mg/kg) prevents cold ischemia-warm reperfusion-induced sinusoidal endothelial cell (SEC) apoptosis and inhibits caspase-3 activation in rat liver by 55 and 94%, respectively.¹ *In vivo*, emricasan reduces alanine aminotransferase (ALT) levels (ED₅₀s = <0.01-0.38 mg/kg) as well as apoptosis and caspase activity in a dose-dependent manner in the α-Fas mouse and D-Gln/LPS rat models of liver injury.² Emricasan reduces caspase-3 and caspase-8 activity, serum ALT levels, hepatocyte apoptosis, and hepatic fibrogenesis in a mouse model of high-fat diet-induced non-alcoholic steatohepatitis (NASH).³ It also enhances islet engraftment and lowers post-transplant fasting glucose levels in a porcine islet autotransplant model.⁴

References

1. Natori, S., Higuchi, H., Contreras, P., *et al.* The caspase inhibitor IDN-6556 prevents caspase activation and apoptosis in sinusoidal endothelial cells during liver preservation injury. *Liver Transpl.* **9(3)**, 278-284 (2003).
2. Hoglen, N.C., Chen, L.S., Fisher, C.D., *et al.* Characterization of IDN-6556 (3-[2-(2-tert-butylphenylaminoxy)amino]-propionylamino]-4-oxo-5-(2,3,5,6-tetrafluoro-phenoxy)-pentanoic acid): A liver-targeted caspase inhibitor. *J. Pharmacol. Exp. Ther.* **309(2)**, 634-640 (2004).
3. Barreyro, F.J., Holod, S., Finocchietto, P.V., *et al.* The pan-caspase inhibitor emricasan (IDN-6556) decreases liver injury and fibrosis in a murine model of non-alcoholic steatohepatitis. *Liver Int.* **35(3)**, 953-966 (2015).
4. McCall, M.D., Maciver, A.M., Kin, T., *et al.* Caspase inhibitor IDN6556 facilitates marginal mass islet engraftment in a porcine islet autotransplant model. *Transplantation* **94(1)**, 30-35 (2012).

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, 09/30/2022

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