

Vorasidenib

Isocitrate dehydrogenase (IDH) 1/2 inhibitor

Vorasidenib is a highly potent, selective, and brain-penetrant inhibitor of mutant isocitrate dehydrogenase 1 and 2 (IDH1/2). It led to >97% inhibition of 2-hydroxyglutarate production in mIDH1 glioma tissue. Vorasidenib significantly improved progression-free survival and delayed time to next intervention in a phase 3 glioma trial.

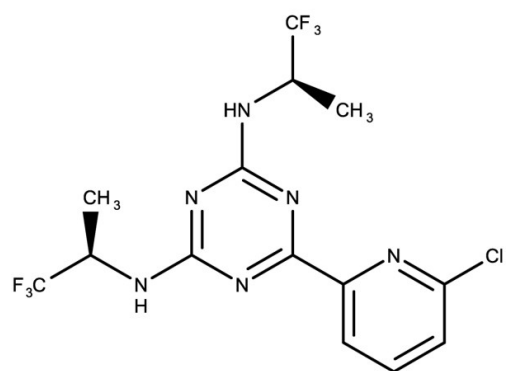
Ordering Information

[Order Online »](#)

ENZ-CHM455-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 6-(6-Chloropyridin-2-yl)-2-N,4-N-bis[(2R)-1,1,1-trifluoropropan-2-yl]-1,3,5-triazine-2,4-diamine

Appearance White solid.

CAS 1644545-52-7

Couple Type Inhibitor

Formula $C_{14}H_{13}ClF_6N_6$

Identity Determined by NMR.

MW 414.74

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (25 mg/ml).

Last modified: January 24, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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