

Venetoclax

Bcl-2 selective inhibitor

Venetoclax is a Bcl-2-selective inhibitor with K_i of 4800-fold more selective versus Bcl-xL and Bcl-w, and no activity to Mcl-1

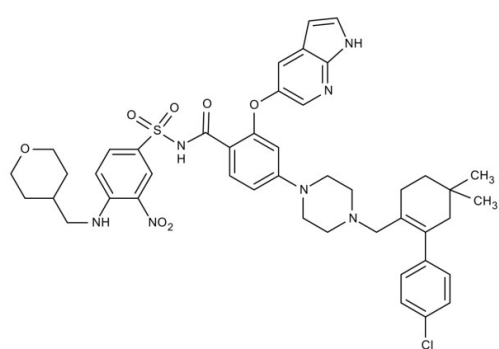
Ordering Information

[Order Online »](#)

ENZ-CHM241-0010	10mg
ENZ-CHM241-0100	100mg
ENZ-CHM241-1000	1g

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Yellow solid.
CAS	1257044-40-8
Couple Target	Bcl-2
Couple Type	Inhibitor
Formula	$C_{45}H_{50}ClN_7O_7S$
MW	868.4
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO.

Last modified: May 29, 2024



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-
be@enzolifesciences.com](mailto:info-be@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)