

Berbamine . dihydrochloride

Berbamine dihydrochloride is an inhibitor of
NF-κB activity

Berbamine dihydrochloride is an inhibitor of NF-κB activity with remarkable anti-myeloma efficacy.

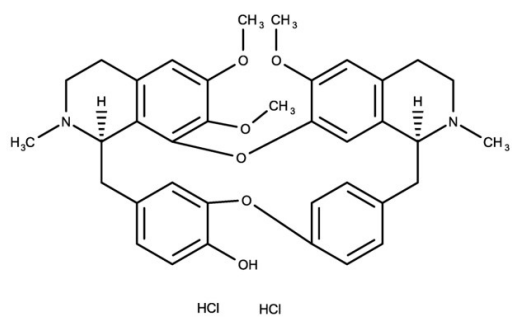
Ordering Information

[Order Online »](#)

ENZ-CHM462-0500	500mg
-----------------	-------

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 16H-1,24:6,9-Dietheno-11,15-metheno-2H-pyrido[2',3':17,18][1,11]dioxacycloeicosino[2,3,4-ij]is oquinolin-12-ol, 3,4,4a,5,16a,17,18,19-octahydro-21,22,26-trimethoxy-4,17-dimethyl-, hydrochl oride (1:2), (4aS,16aR)-

Appearance White to off-white solid.

CAS 6078-17-7

Couple Target NF-kappaB

Couple Type Inhibitor

Formula $C_{37}H_{42}Cl_2N_2O_6$

Identity Determined by NMR.

MW 681.65

Purity ≥96% (HPLC)

Solubility Soluble in DMSO (100mg/ml) or water (25 mg/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-
be@enzolifesciences.com](mailto:info-be@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)