

Bongkreikic acid . triammonium salt

MPTP inhibitor

Bongkreikic acid is an inhibitory ligand of the mitochondrial adenine nucleotide translocase. It inhibits mitochondrial permeability transition pore opening. At 50 μ M it inhibits dexamethasone- apoptosis in thymocytes, preventing the PARP cleavage and DEVDase (caspase-3 like) activity normally produced by these treatments.

Citations: 3

[View Online »](#)

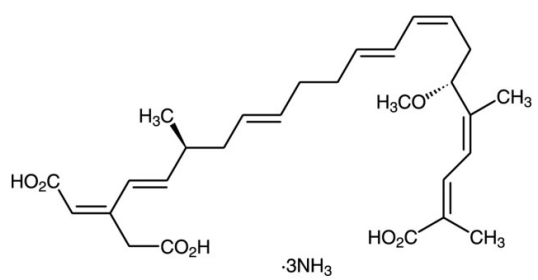
Ordering Information

[Order Online »](#)

BML-CM113-1000	1mg
----------------	-----

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

Appearance Beige to pink lyophilized solid.

CAS 1177154-51-6

Couple Target MPTP

Couple Type Inhibitor

Formula $C_{28}H_{38}O_7 \cdot 3 NH_3$

MW 537.7

Purity ≥92% (HPLC)

Reconstitution Reconstitute to 1mg/ml with distilled water

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)