

BML-277

Chk inhibitor

BML-277 is a highly selective inhibitor of Chk2 ($IC_{50}=15nM$). It displays potent radioprotective activity and prevents apoptosis of human T cells subjected to ionizing radiation ($EC_{50}=3-7.6\mu M$). Useful tool for dissecting the role of Chk2 in cellular signaling.

Citations: 3

[View Online »](#)

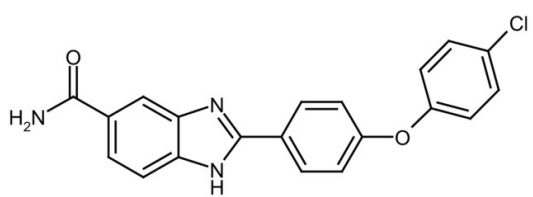
Ordering Information

[Order Online »](#)

BML-EI388-0010	10mg
----------------	------

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

Alternative Name Chk2 Inhibitor II

Appearance Off-white solid.

CAS 516480-79-8

Couple Target Chk

Couple Type Inhibitor

Formula $C_{20}H_{14}ClN_3O_2$

MW 363.8

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (20mg/ml) or ethanol (20mg/ml).

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-
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)