

BCTC

TRPV1 antagonist

BCTC is a potent and selective inhibitor of capsaicin-induced ($IC_{50}=35$ nM) and acid-induced ($IC_{50}=6$ nM) VR_1 activation. It displays analgesic activity in rat models of inflammatory and neuropathic pain. TRPV1 antagonist.

Citations: 2

[View Online »](#)

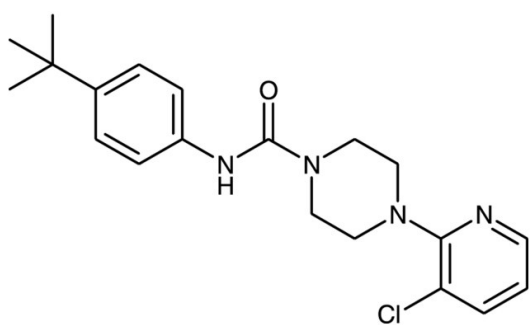
Ordering Information

[Order Online »](#)

BML-CA231-0010	10mg
BML-CA231-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Tan solid

CAS 393514-24-4

Couple Target TRP channel

Couple Type Inhibitor

Formula $C_{20}H_{25}ClN_4O$

MW 372.9

MeltingPoint 150-152°C

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (20 mg/ml) or ethanol (15 mg/ml with warming).

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)