

Butyrolactone I

CDK inhibitor

Cell permeable and selective inhibitor of the cyclin-dependent kinases CDK1/cyclin B, CDK2 and CDK5. Inhibits cell cycle progression at the G1/S and G2/M transitions. G1/S and G2/M transitions, in W138 cells, were inhibited, concurrent with inhibition of the phosphorylations of, respectively, retinoblastoma protein and histone H1. It is an important tool for probing the cellular roles of CDKs. Blocks Fas-induced apoptosis. Displays antitumor activity.

Citations: 15

[View Online »](#)

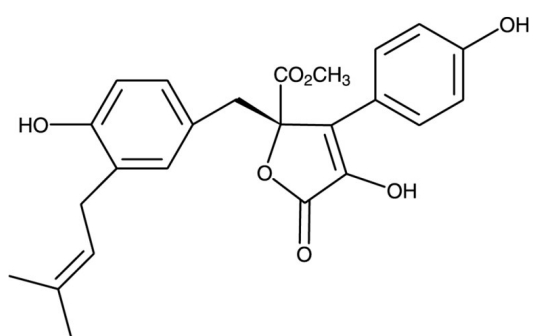
Ordering Information

[Order Online »](#)

BML-CC210-1000	1mg
BML-CC210-0200	200µg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	87414-49-1
Couple Target	CDK
Couple Type	Inhibitor
Formula	$C_{24}H_{24}O_7$
MW	424.5
Purity	≥99% (TLC)
Solubility	Soluble in 100% ethanol (25mg/ml) or DMSO(>10mg/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-
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)