

5,6-Dichloro-1- β -D-ribofuranosylbenzimidazole

Casein kinase inhibitor

DRB is a potent and specific inhibitor of casein kinase II (CKII, $IC_{50} \sim 6 \mu M$). It has been used to inhibit RNA polymerase II mediated-transcription which may be dependent on CKII and its interaction with ATF-1.

Citations: 10

[View Online »](#)

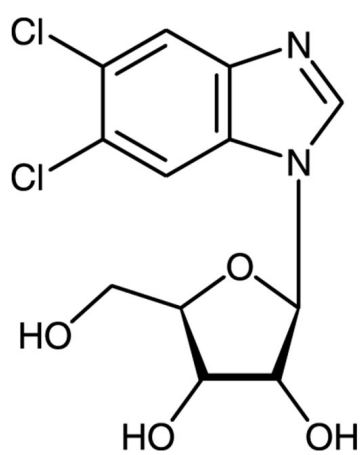
Ordering Information

[Order Online »](#)

BML-EI231-0050	50mg
----------------	------

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.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DRB, 5,6-Dichlorobenzimidazole riboside
Appearance	White solid.
CAS	53-85-0
Couple Target	Casein kinase
Couple Type	Inhibitor
Formula	$C_{12}H_{12}Cl_2N_2O_4$
Identity	Determined by 1H-NMR.
MW	319.1
Purity	≥97% (TLC)
Solubility	Soluble in DMSO (20mg/ml), 100% ethanol, and dimethyl formamide (100mM).

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)