

DRB

Inhibitor of Casein kinase II

Inhibitor of casein kinase II and of RNA synthesis in eucaryotic cells.

BLG-D012-500 (5 x 100 µmol pack size) is not sold in the U.S. or Canada.

Please [contact us](#) for available options.

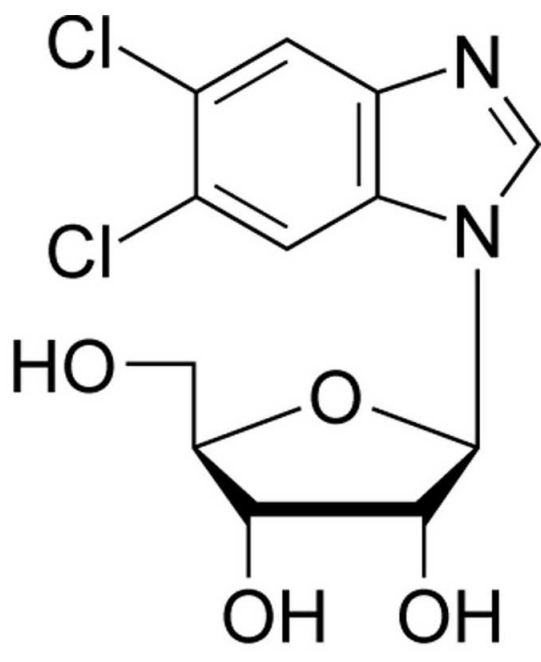
Ordering Information

[Order Online »](#)

BLG-D012-500	5x100µmol
--------------	-----------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5,6-Dichloro-1-β-D-ribofuranosyl benzimidazole

CAS 53-85-0

Couple Type Modified nucleotides

Formula $C_{12}H_{12}Cl_2N_2O_4$

MW 319.1

Purity > 98% HPLC

Quantity 100μmol≈32mg

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com