

# CDK1/Cyclin B (human), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

|                |      |
|----------------|------|
| BML-SE295-0010 | 10µg |
|----------------|------|

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Cyclin-dependent kinase 1/Cyclin B, p34cdc2/Cyclin B

**Application Notes** Useful for kinetic and functional studies, phosphorylation of target substrates, drug screening.

**Formulation** Liquid. In 50 mM HEPES, pH 7.5, containing 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, and 20% glycerol.

**Source** Produced in Baculovirus infected Sf9 insect cells.

**UniProt ID** P06493 (CDK1)

**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)