

# CDK2/Cyclin E (human), (recombinant) (GST- tag)

Active serine/threonine kinase, using histone H1 as substrate.

**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 2/Cyclin B

**Application Notes**

Useful for studying enzyme kinetics, phosphorylation of target substrates, and drug screening.

**Formulation**

Liquid. In 50mM TRIS-HCl, pH 8.0, containing 100mM sodium chloride, 5mM DTT, 15mM reduced glutathione and 20% glycerol.

**MW**

60.2 kDa (CDK2) 72.0 kDa (cyclin E)

**Source**

Produced in insect cells. GST fusion proteins containing CDK2 (aa 1-298) and residues cyclin E (aa 1-394).  
Produced in a baculovirus expression system.

**UniProt ID**

P24941 (CDK2)

**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)