

# CDK5/p25 (human), (recombinant) (GST- tag)

CDK5 is the most abundant member of the cyclin-dependent kinase family in the mammalian brain. It exists as a heterodimer of CDK5 and p25, which is necessary for kinase activity. It is a crucial regulator of neuronal migration in the developing CNS.

## Ordering Information

[Order Online »](#)

BML-SE471-0005	5µg
----------------	-----

## 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 5

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

MW      59 and 49 kDa

Source      Produced in insect cells. Full length active enzyme. Produced in a baculovirus expression system.

UniProt ID      Q00535



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)