

# CDC25C (human), (recombinant)

CDC25 protein tyrosine phosphatases activate cyclin dependent kinases by dephosphorylating them, promoting cell division. All three of the genes encoding the CDC25 phosphatases are considered potential oncogenes and the enzymes are drug discovery targets. CDC25C activates CDC2/cyclin B which controls entry into mitosis. CDC25C, itself, is regulated by phosphorylation and binding to 14-3-3 proteins.

## Ordering Information

[Order Online »](#)

<b>BML-SE366-0050</b>	50µg
-----------------------	------

## 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**      Cell division cycle 25C, Dual specificity phosphatase CDC25C

**Application Notes**      Useful for kinetic studies and high throughput screening.

**Formulation**      Liquid. In 50mM TRIS/HCl, pH 8.0, 150mM NaCl, 5mM DTT, 0.1mM EDTA, 0.03% Brij 35, 20% glycerol.

**MW**      54 kDa

**Purity Detail**      Partially purified by single-step affinity chromatography and gel filtration.

**Source**      Produced in *E. coli*. Full length human CDC25C.

**Specific Activity**      ≥14 U/mg assayed by 3-O-methyl fluorescein phosphate (OMFP) hydrolysis at pH 8.2, 30°C. One unit is equal to 1 nmole phosphate hydrolyzed from OMFP per minute.

**UniProt ID**      P30307



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)