

CDC25B (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. In mammalian cells, the CDC25B isoform activates cyclin A- and cyclin B1-containing complexes and is necessary for entry into mitosis. CDC25B ^{-/-} female mice are sterile because their oocytes are in permanent meiotic arrest resulting from the inability to activate CDK11. CDC25B shuttles between the nucleus and the cytoplasm. The protein is in the nucleus during M and G1 and moves to the cytoplasm during S and G2.

Ordering Information

[Order Online »](#)

| | |
|-----------------------|------|
| BML-SE365-0050 | 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 25B, Dual specificity phosphatase CDC25B

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.

Gene/Protein Identifier NM_021873 (RefSeq)

MW 65 kDa

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

Source Produced in *E. coli*. Full length human CDC25B.

Specific Activity ≥10 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 P30305



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)