

# PDE4B2 (catalytic domain) (human), (recombinant)

PDE4B interacts directly with DISC1, which is a major risk factor associated with schizophrenia<sup>1</sup>. PDE4B hydrolyzes cAMP and DISC1 may regulate this activity.

## Ordering Information

[Order Online »](#)

|                       |      |
|-----------------------|------|
| <b>BML-SE522-0020</b> | 20µg |
|-----------------------|------|

## 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

Phosphodiesterase 4B2

Application Notes

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

MW

37 kDa

Source

Produced in *E. coli*. Active catalytic domain of phosphodiesterase 4B2.

UniProt ID

Q07343#Q07343-2

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)