

# PDE7A (catalytic domain) (human), (recombinant)

PDE7A is expressed in immune and pro-inflammatory cells. It is being investigated as an alternate target to PDE4 for the inhibition of inflammation that is associated with many inflammatory diseases including asthma, chronic obstructive pulmonary disease, atopic dermatitis, psoriasis, lupus, rheumatoid arthritis and multiple sclerosis.

## Ordering Information

[Order Online »](#)

BML-SE524-0020	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 7A

MW      39 kDa

Source      Produced in *E. coli*. Active catalytic domain of phosphodiesterase 7A.

UniProt ID      Q13946



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)