

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.

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 <i>E. coli</i> . Active catalytic domain of phosphodiesterase 7A.
UniProt ID	Q13946

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com