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

Produced in E. coli. Active catalytic domain of Source

phosphodiesterase 7A.

UniProt ID Q13946

Last modified: May 29, 2024

info-