

PDE10A (catalytic domain) (human), (recombinant)

This enzyme can hydrolyze both cAMP and cGMP, but has a higher affinity for cAMP. In the brain it is present in neuronal cell bodies and neuropil of the striatum, in striatonigral and striatopallidal white matter tracks, and in the substantia nigra and globus pallidus. Outside the brain it is found in the testis, epididymal sperm, and enteric ganglia¹. It is under investigation as a drug target for schizophrenia and diabetes.

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 10A
Application Notes	Useful for kinetic and functional studies, phosphorylation of target substrates, drug screening.
Formulation	Liquid. In 20mM Tris-HCl, pH 7.5, containing 150mM NaCl, 2mM DTT.
MW	37 kDa
Source	Produced in <i>E. coli</i> . Active catalytic domain of phosphodiesterase 10A.
UniProt ID	Q9Y233

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