

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.

Ordering Information

[Order Online »](#)

BML-SE525-0020	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 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 *E. coli*. 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](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)