

PDE5A (human), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

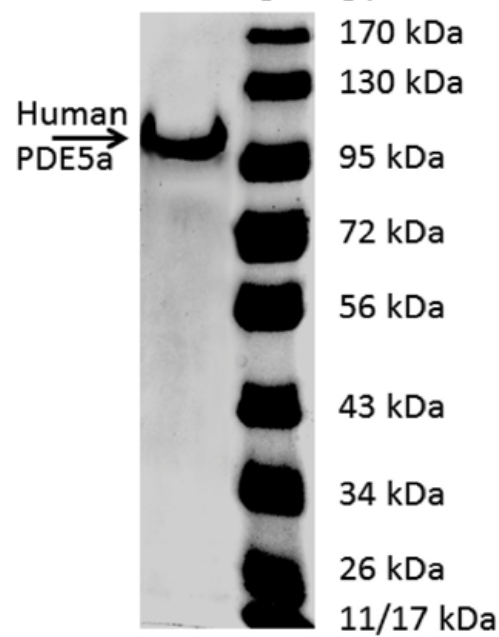
[Order Online »](#)

ENZ-PRT126-0010	10µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Coomassie staining, 5 ug purified hPDE5A



Handling & Storage

Use/Stability Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Phosphodiesterase 5A, cGMP-specific 3',5'-cyclic phosphodiesterase, CGB-PDE

Formulation Liquid. In 50mM triethanolamine, pH 7.4, containing 1mM DTT, 100mM NaCl, 20% glycerol, 0.5mM EDTA.

Purity ≥95%

Source Produced in Sf9 insect cells. Recombinant human PDE5A (phosphodiesterase 5A) is fused at the C-terminus to a His-tag. Produced in a baculovirus expression system.

UniProt ID O76074

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)