

Pregnancy-associated plasma protein A (human) monoclonal antibody (PAPP A 1-1)

Pregnancy-associated plasma protein-A (PAPP-A, Insulin-like growth factor binding protein-4 protease, PAPP-A-proMBP complex) is a large heterotetrameric glycoprotein of approximately 500kDa, which was first discovered in serum from pregnant women. The heterotetramer consists of two PAPP-A subunits and two proMBP subunits. Low maternal serum levels of PAPP-A in first trimester biochemical screening is used as a marker of Down's syndrome (trisomy 21).

This antibody is covered by our [Worry-Free Guarantee](#).

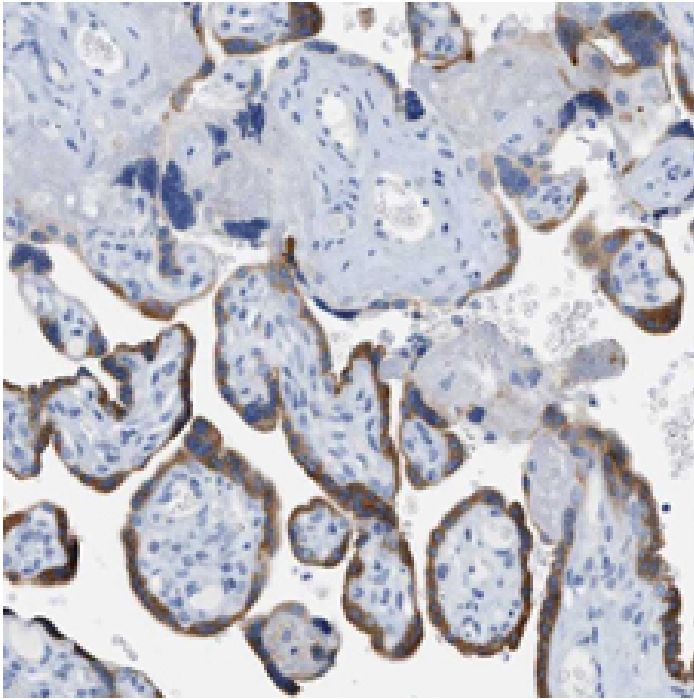
Ordering Information

[Order Online »](#)

BPD-ABS-006-01-1	1mg
BPD-ABS-006-01-04	400µg

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical analysis of placental tissue using PAPP-A (human), mAb (1_1).

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name PAPP-A, Pappalysin-1, Insulin-like growth factor binding protein-4 protease

Application ELISA, IHC, WB

Clone PAPP A 1-1

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

GenBank ID 5069

Host Mouse

Immunogen Purified human PAPP-A.

Isotype IgG1κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human

Specificity Specific for native human PAPP-A.

UniProt ID Q13219

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



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-
fr@enzolifesciences.com](mailto:info-fr@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)