

Pregnancy-associated plasma protein A (human) monoclonal antibody (9)

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 hetreotetramer 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-BTE-004-09-1	1mg
BPD-BTE-004-09-04	400µg

Manuals, SDS & CofA

[View Online »](#)

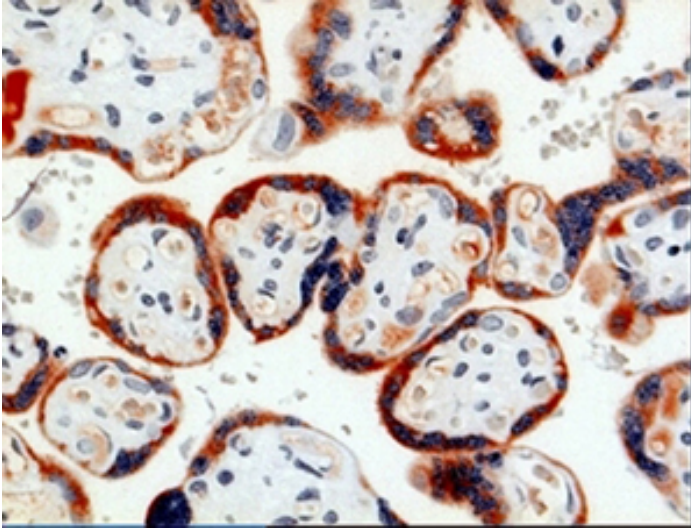


Figure 2: Immunohistochemical analysis of human placenta tissue using PAPP-A (human), mAb specific for human pregnancy associated plasma protein A (PAPP-A).

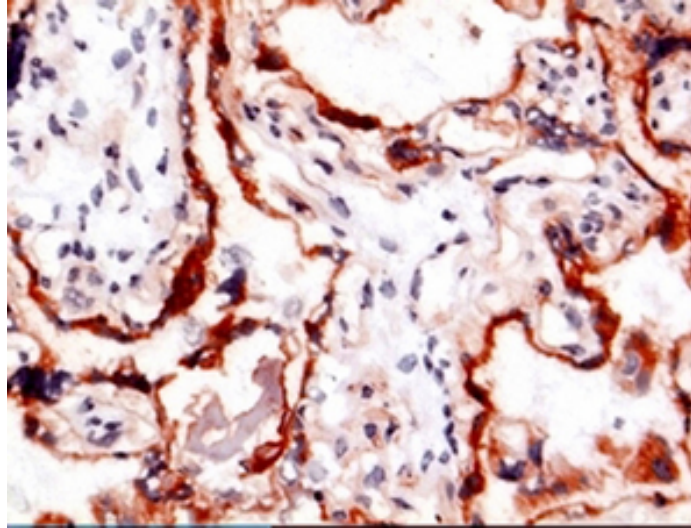


Figure 1: Immunohistochemical analysis of human placenta tissue using PAPP-A (human), mAb specific for human pregnancy associated plasma protein A (PAPP-A).

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 (FS), IHC (PS), WB
Clone	9
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	IgG2bk
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Human
Specificity	Specific for native human PAPP-A. Heat and SDS denatured PAPP-A is also recognized, but not acid treated 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-
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)