

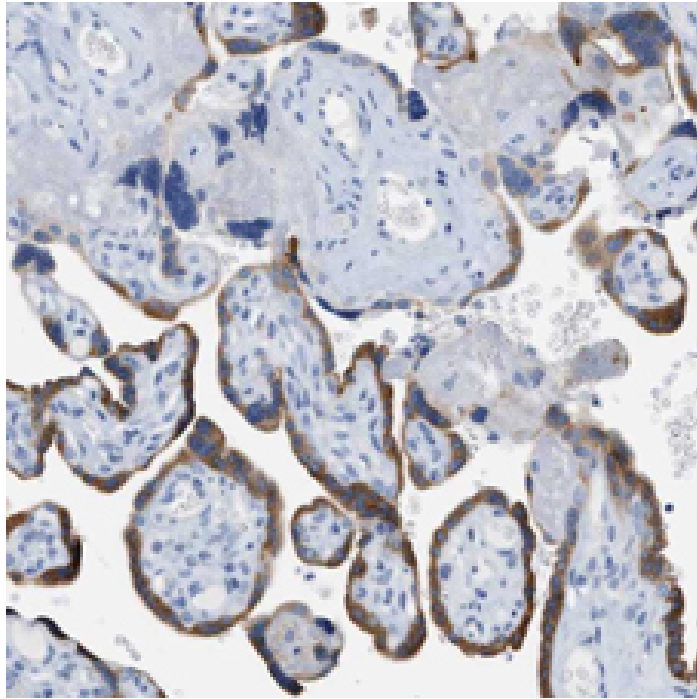
Pregnancy-associated plasma protein A (human) monoclonal antibody (PAPP A 1-1)

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](#).

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

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

Last modified: May 29, 2024