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

cy-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-ABS-006-24-1	1mg
BPD-ABS-006-24-04	400μg

Manuals, SDS & CofA

**View Online** »

## **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 1-24

**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 IgG1ĸ

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com