

Anti-PLAC3

Rabbit Polyclonal Antibody

Catalog # P38-363CR

Lot # O2361-43

Cited Applications

E, WB, ICC, IF

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the PLAC3 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

PLAC3 antibody was raised against a 15 amino acid synthetic peptide near the amino terminus of human PLAC3

Formulation

PBS + 0.02% sodium azide.

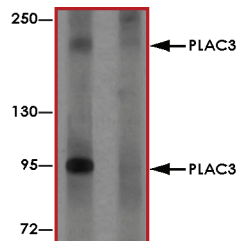
Stability

1yr at -20°C from date of shipment.

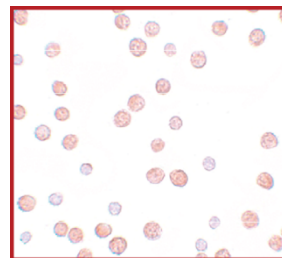
References

1. Overgaard MT et al: Pregnancy-associated plasma protein-A2 (PAPP-A2), a novel insulin-like growth factor-binding protein-5 proteinase. *J. Biol. Chem.*2001; 276:1849-53.
2. Wang J et al: Expression of pregnancy-associated plasma protein A2 during pregnancy in human and mouse. *J. Endocrin.*2009; 202:337-45.
3. Nishizawa H et al: Increased levels of pregnancy-associated plasma protein-A2 in the serum of pre-eclamptic patients. *Mol. Hum. Reprod.*2008; 14:595-602.
4. Buimer M et al: Seven placental transcripts characterize HELLP syndrome. *Placenta*2008; 29:444-53.

Sample Data



Western blot analysis of PLAC3 in HeLa cell lysate with PLAC3 antibody at 1 µg/mL in (A) the absence and (B) the presence of blocking peptide.



Immunocytochemistry of PLAC3 in HeLa cells with PLAC3 antibody at 10 µg/mL.

Anti-PLAC3

Rabbit Polyclonal Antibody

Catalog Number

P38-363CR

Specific Lot Number

O2361-43

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.