APC-AAT complexes, neoepitope (human) monoclonal antibody (PC 7)

Protein C is a vitamin K-dependent serine protease produced in the liver and made up of 2 polypeptide chains. The 62kDa proenzyme is activated by thrombin, and the active enzyme cleaves factor Va and VIIIa, thus inhibiting blood coagulation. The molecular weight of the active enzyme is 55kDa. The normal concentration in human plasma is approximately 1-3 ng/mL due to the very fast turnover and the proenzyme concentration of approximately 3 ug/mL. The activated protein C (APC) is inhibited by members of the serine protease inhibitor (serpin) family, of which alpha1-antitrypsin (AAT) and protein C inhibitor (PCI) are the most important.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BPD-ABS-001-07-1	1mg
BPD-ABS-001-07-04	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 Activated protein C and α1-antitrypsin complex

Application ELISA

Clone PC 7

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M

sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Recombinant human activated protein C.

lsotype lgG1κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human

Specificity Specific for a conformation-dependent neoepitope that is

expressed in activated protein C upon complex-formation with α 1-antitrypsin. No reaction is seen to non-complexed α 1-antitrypsin and only very little cross reaction to protein C zymogen. Note that specificity is calcium dependent.

UniProt ID P04070 (APC), P01009 (AAT)

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com