

# 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](#).

**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 $\alpha$ 1-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.
Isotype	IgG1 $\kappa$
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 $\alpha$ 1-antitrypsin. No reaction is seen to non-complexed $\alpha$ 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 <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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)