

Anti-Amyloid beta

Rabbit Polyclonal Antibody

Catalog # **A06-563R**

Lot # Z2014-1

Cited Applications

WB, ELISA, IHC, IF

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

Specificity

Recognizes the amyloid beta protein

Cross Reactivity

Human and Mouse

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Amyloid beta antibody was raised against a synthetic peptide corresponding to the amino terminus (aa 1-14) of human beta amyloid conjugated to KLH using maleimide.

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) sodium azide

Stability

1yr at -20°C from date of shipment

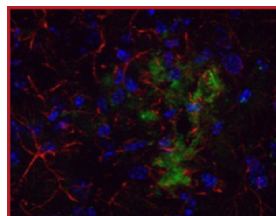
Scientific Background

Amyloid beta, often abbreviated as A-beta, is a protein that builds up in the brains of persons with Alzheimer's disease, collecting in clumps called plaques or senile plaques. While some researchers question whether beta amyloid is the cause of the dementia, most agree that it is involved in the disruption of thinking that is a hallmark of the disease. In some cases of familial Alzheimer's disease, mutations in genes for the proteins called the presenilins lead to increased production of amyloid. Researchers have been looking at how presenilin-1 in particular contributes to the excess buildup of beta amyloid. Presenilin-1 apparently acts to increase the activity of gamma-secretase, an enzyme that changes a normal protein (amyloid precursor protein or APP) into beta amyloid itself. Furthermore, presenilin-1 might be gamma-secretase.

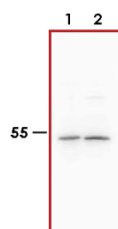
References

1. Frenkel D, N-terminal EFRH sequence of Alzheimer's beta-amyloid peptide represents the epitope of its anti-aggregating antibodies.; J Neuroimmunol, 1998 Aug 1
2. Bonaiuto C , Activation of nuclear factor-kappa B by beta-amyloid peptides and interferon-gamma in murine microglia; J Neuroimmunol, 1997 Jul
3. Suo Z, Alzheimer's beta-amyloid peptides induce inflammatory cascade in human vascular cells: the roles of cytokines and CD40; Brain Res, 1998 Oct 5

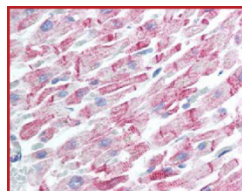
Sample Data



Immunohistochemical detection of Amyloid beta using Anti-Amyloid beta (1:200) on TG APP23 mouse brain cortex frozen sections.



Mouse Brain (Lane 1) and Mouse Spinal Cord (Lane 2) were run on a 4-20% gradient gel, and probed with Anti-Amyloid beta 1:1000



Human Heart (formalin-fixed, paraffin-embedded) stained with Anti-Amyloid beta Antibody at 5 µg/ml

Anti-Amyloid beta

Rabbit Polyclonal Antibody

Catalog #

A06-563R

Lot #

Z2014-1

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.