

Ub+1 Polyclonal Antibody

(Catalog No. A-8504)

Background

Ub+1 is a novel form of ubiquitin that has been selectively observed in the brains of Alzheimer's patients, including those with nonfamilial Alzheimer's disease. Ub⁺ is an efficient substrate for polyubiquitinylation *in vitro* and in transfected human cells. The resulting polyubiquitin chains are refractory to disassembly by deubiquitinating enzymes and potently inhibit the degradation of a polyubiquitinated substrate by purified 26S proteasomes. Thus, expression of Ub⁺ in aging brain could result in dominant inhibition of the Ub-proteasome system, leading to neuropathologic consequences

Description

Rabbit polyclonal antibody to Ub+1.

Formulation

Liquid. Antiserum containing 0.1% sodium azide.

Immunogen

Synthetic peptide corresponding to aa 76-95 of human Ub+1.

Specificity

Human

Purification

Serum

Storage

Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

Ubiquitin+1

Application

IHC, WB (1:1000)

Ordering Information

Products	Size	Cat. No.
Ub+1 Polyclonal Antibody	100 µl	A-8504-100



▲ Western blot analysis of Ub+1 (1 µg) after SDS- PAGE followed by blotting onto PVDF and probing with Ub+1 Polyclonal Antibody with [lane (a)] and without [lane (b)] peptide blockade. Antibody dilution 1:1000 using ECL procedure

This product is for research purposes only. Not intended for use in diagnostic procedures.