

---

# ATG4A Polyclonal Antibody, FITC Conjugated

(Catalog # A51135)

---

## Background

Cysteine protease required for the cytoplasm to vacuole transport (Cvt) and autophagy. Cleaves the C-terminal amino acid of ATG8 family proteins to reveal a C-terminal glycine. Exposure of the glycine at the C-terminus is essential for ATG8 proteins conjugation to phosphatidylethanolamine (PE) and insertion to membranes, which is necessary for autophagy. Preferred substrate is GABARAPL2 followed by MAP1LC3A and GABARAP. Has also an activity of delipidating enzyme for the PE-conjugated forms.

## Description

ATG4A Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

## Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

## Specificity

Human

## Isotype

IgG

## Uniprot ID

Q8WYN0

## Purification

>95%, Protein G purified

## Immunogen

Recombinant Human Cysteine protease ATG4A protein (1-214AA)

## Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

## Alternative Names

AUT-like 2 cysteine endopeptidase, Autophagin-2, Autophagy-related cysteine endopeptidase 2, Autophagy-related protein 4 homolog A, hAPG4A, ATG4A, APG4A, AUTL2

## Application

ELISA; Recommended dilution: ELISA 1:100-1:500

---