

PRODUCT INFORMATION



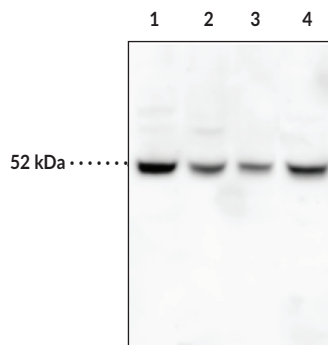
Beclin 1 Monoclonal Antibody (Clone 5F7)

Item No. 25964

Overview and Properties

Contents: This vial contains 100 µg of protein G-purified monoclonal antibody.
Synonyms: BECN1, Coiled-coil Myosin-like Bcl-2-interacting Protein, Protein GT197
Immunogen: Full-length recombinant human Beclin 1
Species Reactivity: (+) Human and murine; other species not tested
Uniprot No.: Q14457
Form: Liquid
Storage: -20°C (as supplied)
Stability: ≥3 years
Storage Buffer: PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Clone: 5F7
Host: Mouse
Isotype: IgG2b
Applications: ELISA and Western blot (WB); the recommended starting dilution for ELISA and WB is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: RAW 264.7 Lysate (50 µg)
Lane 2: HepG2 Lysate (50 µg)
Lane 3: HEK293 Lysate (50 µg)
Lane 4: A431 Lysate (50 µg)

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/06/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Beclin 1 (BECN1) is a core component of the class III phosphatidylinositol 3-kinase (PI3K) complex, which generates phosphatidylinositol 3-phosphate (PI3P) and has functions in macroautophagy and the endocytic pathway when associated with ATG14 and UVRAG, respectively.¹ BECN1 is comprised of a BH3 domain involved in Bcl-2/Bcl-xL binding, a coiled-coil domain, and an evolutionarily conserved domain (ECD) that is required for association with the PI3K complex.^{2,3} Binding of BECN1 to various BECN1-interacting proteins can regulate the activity of the ATG14-containing PI3K complex, which is required for nucleation of the phagophore during macroautophagy.^{1,4} Binding of the anti-apoptotic protein Bcl-2 to BECN1 prevents the association of BECN1 with the PI3K complex and inhibits autophagy, whereas AMBRA1, which also binds directly to BECN1, positively regulates the PI3K complex to facilitate autophagy.^{1,3} Transfection of autophagy-deficient MCF-7 human breast cancer cells with wild-type BECN1, but not a BECN1 mutant lacking the ECD, increases starvation-induced autophagy *in vitro* and inhibits tumor growth in an MCF-7 mouse xenograft model, indicating that BECN1 has ECD-dependent tumor suppressor activity.² BECN1 heterozygous mice expressing human amyloid precursor protein (*APP⁺Becn1^{+/-}* mice) have decreased autophagy in the cerebral cortex, increased extracellular amyloid- β (A β) deposits in the frontal cortex, and increased intraneuronal A β accumulation in the hippocampus and frontoparietal cortex compared to *APP⁺Becn1^{+/+}* mice.⁵ BECN1 may also have roles in ischemia/reperfusion injury, Niemann-Pick type C disease, and protection against viral infection.³ Cayman's Beclin 1 Monoclonal Antibody (Clone 5F7) can be used for Western blot and ELISA applications. The antibody recognizes Beclin 1 at approximately 52 kDa from human samples.

References

1. Parzych, K.R. and Klionsky, D.J. An overview of autophagy: Morphology, mechanism, and regulation. *Antioxid. Redox Signal.* **20(3)**, 460-473 (2014).
2. Furuya, N., Yu, J., Byfield, M., *et al.* The evolutionarily conserved domain of Beclin 1 is required for Vps34 binding, autophagy and tumor suppressor function. *Autophagy* **1(1)**, 46-52 (2005).
3. Cao, Y. and Klionsky, D.J. Physiological functions of Atg6/Beclin 1: A unique autophagy-related protein. *Cell Res.* **17(10)**, 839-849 (2007).
4. Burman, C. and Ktistakis, N.T. Regulation of autophagy by phosphatidylinositol 3-phosphate. *FEBS Lett.* **584(7)**, 1302-1312 (2010).
5. Pickford, D., Masliah, E., Britschgi, M., *et al.* The autophagy-related protein beclin 1 shows reduced expression in early Alzheimer disease and regulates amyloid β accumulation in mice. *J. Clin. Invest.* **118(6)**, 2190-2199 (2008).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM