

PRODUCT INFORMATION



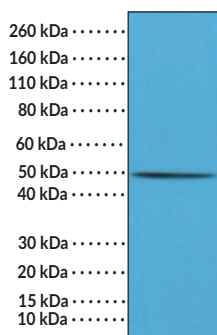
BAG-1L Rabbit Monoclonal Antibody (Clone RM310)

Item No. 32261

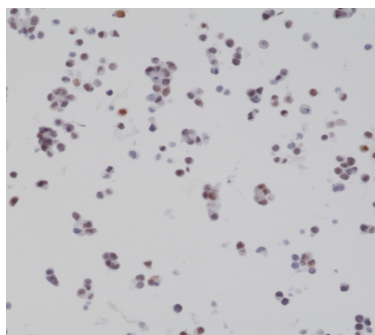
Overview and Properties

Contents:	This vial contains 100 μ l of protein A-affinity purified monoclonal antibody.
Synonyms:	BAG Co-chaperone 1L, BAG Family Molecular Chaperone Regulator 1L, Bcl-2-associated Anthanogene-1L
Immunogen:	Peptide corresponding to a region near the N-terminus of human BAG-1L
Cross Reactivity:	(+) BAG-1L; (-) BAG-1M, BAG-1S
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM310
Host:	Rabbit
Isotype:	IgG
Applications:	Immunocytochemistry (ICC), Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:200-1:1,000 for ICC, 1:500-1:2,000 for IHC, and 1:1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

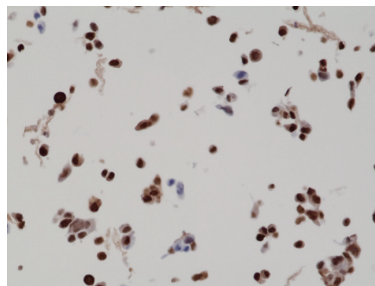
Images



WB of HeLa cell lysate using BAG-1L Rabbit Monoclonal Antibody (Clone RM310) at a 1:1,000 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded 22RV1 cells using BAG-1L Rabbit Monoclonal Antibody (Clone RM310) at a 1:2,000 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded BAG-1L overexpressing LNCaP cells using BAG-1L Rabbit Monoclonal Antibody (Clone RM310) at a 1:2,000 dilution.

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, 01/30/2024

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

Bcl-2-associated anthanogene 1 (BAG-1) is a multifunctional co-chaperone protein with roles in apoptosis and transcription, as well as cell motility, survival, and proliferation.^{1,2} It is encoded by a single gene, *BAG1*, that produces several isoforms with variable N-termini through alternate translation initiation of a single mRNA transcript.³ BAG-1L, the longest isoform, is localized to the nucleus and composed of a nuclear localization sequence, a basic DNA-binding domain, a ubiquitin-like domain, and an Hsp70- and Hsc70-binding domain.¹ It functions as a co-chaperone to enhance transactivation of the androgen receptor *via* interaction with the partially folded and intrinsically disordered N-terminal tau-5 domain, which stabilizes the receptor to facilitate chromatin binding and activates gene transcription.⁴ Overexpression of BAG-1L inhibits apoptosis induced by cisplatin (Item No. 13119) in HeLa cells stably transfected with pEZ-M02-BAG-1L plasmid DNA, and BAG-1L levels are increased in tumors isolated from patients with castration-resistant prostate cancer (CRPC).^{3,4} Cayman's BAG-1L Rabbit Monoclonal Antibody (Clone RM310) can be used for immunocytochemistry (ICC), immunohistochemistry (IHC), and Western blot (WB) applications.

References

1. Gehring, U. Activities of the cochaperones Hap46/BAG-1M and Hap50/BAG-1L and isoforms. *Cell Stress Chaperones* **11**(4), 295-303 (2006).
2. Dobbyn, H.C., Hill, K., Hamilton, T.L., *et al.* Regulation of BAG-1 IRES-mediated translation following chemotoxic stress. *Oncogene* **27**(8), 1167-1174 (2008).
3. Ozfiliz, P., Arisan, E.D., Coker-Gurkan, A., *et al.* Bag-1L is a stress-withstand molecule prevents the downregulation of Mcl-1 and c-Raf under control of heat shock proteins in cisplatin treated HeLa cervix cancer cells. *Asian Pac. J. Cancer Prev.* **15**(11), 4475-4482 (2014).
4. Cato, L., Neeb, A., Sharp, A., *et al.* Development of Bag-1L as a therapeutic target in androgen receptor-dependent prostate cancer. *Elife* **6**, e27159 (2017).

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