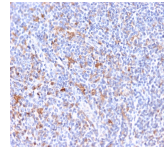


RCAS1 / Estrogen Receptor Binding Site Associated, Antigen 9 Monoclonal Mouse Antibody (CPTC-EBAG9-2)



Product Description

EBAG9, also known as RCAS1, is an estrogen-transcribed protein. Soluble and membranous RCAS1 proteins may play a role in the immune escape of tumor cells by promoting T lymphocyte inhibition of growth and apoptosis. RCAS1 is expressed in a wide variety of cancers, including uterine, ovarian, and lung cancer cells, and acts as a ligand for a putative receptor present on peripheral lymphocytes. RCAS1 is highly expressed not only in cancer cells but also in non-tumor bile duct cells subject to immune attack. RCAS1 inhibits the in vitro growth of receptor-expressing cells and induces apoptosis, contributing to the ability of tumor cells to evade host immune surveillance. High expression of RCAS1 significantly correlates with tumor progression and with poor outcome for many cancer patients.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Product attributes

Antibody number	#2502
Antibody reactivity (target)	Estrogen Receptor Binding Site Associated, Antigen 9, RCAS1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-EBAG9-2
Isotype	IgG2a, kappa
Molecular weight	32 kDa
Synonyms	BAG9; cancer associated surface antigen; cancer associated surface antigen RCAS1; EB9; estrogen receptor binding fragment associated gene 9; PDAF; receptor binding cancer antigen expressed on SiSo cells
Human gene symbol	EBAG9
Entrez gene ID	9166
SwissProt	O00559
Unigene	409368
Immunogen	Recombinant full-length human EBAG9 protein
Verified antibody applications	IHC (FPPE) (verified), WB (verified)
Antibody target cellular localization	Golgi apparatus
Species reactivity	Human
Positive control	MCF-7, Jurkat; 293 cell lysates; Human skeletal muscle tissue.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and of LI-COR Bioscience.

This datasheet was generated on September 1, 2024 at 08:25:11 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/rcas1-estrogen-receptor-binding-site-associated-antigen-9-monoclonal-mouse-antibody-cptc-ebag9-2/>