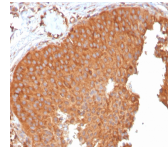


# PAI-RBP1 / SERBP1 / SERPINE1 Monoclonal Mouse Antibody (SERBP1/3496)



## Product Description

PAI-RBP1 (plasminogen activator inhibitor 1 RNA-binding protein), also known as SERBP1 (SERPINE1 mRNA-binding protein 1), CGI-55, CHD3IP (chromodomain helicase DNA binding protein 3 interacting protein), HABP4L or PAI-RBP1, is a membrane-associated protein that localizes to the nucleus, the perinuclear region of the cytoplasm and the plasma membrane. PAI-RBP1 is believed to play a role in the regulation of mRNA stability, as it specifically binds to the CRS (cyclic nucleotide-responsive sequence) motif of the PAI-1 mRNA and acts to stabilize the mRNA and regulate its expression. In addition, PAI-RBP1 interacts with Mi2- and may be involved in chromatin remodeling. PAI-RBP1 also interacts with PGRMC1 and participates in the transduction of Progesterone s antiapoptotic action in ovarian cell types. The gene encoding PAI-RBP1 is overexpressed in ovarian cancer, suggesting a possible role for PAI-RBP1 in tumorigenesis and tumor metastasis. 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

## Product attributes

Antibody number	#3496
Antibody reactivity (target)	PAI-RBP1, SERBP1, SERPINE1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	SERBP1/3496
Isotype	IgG2c, kappa
Molecular weight	60 kDa
Synonyms	CGI-55; CHD3IP; Chromodomain helicase DNA binding protein 3 interacting protein; HABP4L; PAI1 RNA-binding protein 1; PAIRBP1; Plasminogen activator inhibitor 1 RNA binding protein; SERPINE1 mRNA binding protein 1
Human gene symbol	SERBP1
Entrez gene ID	26135
SwissProt	Q8NC51
Unigene	530412
Immunogen	Recombinant fragment of human SERBP1 protein (around aa3-139) (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Nucleus & cytoplasm, Perinuclear
Species reactivity	Human
Positive control	Human kidney or bladder tissue (IHC).
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
Validated in protein array	Monospecific
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	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
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, LI-COR Bioscience.

This datasheet was generated on June 29, 2024 at 09:20:35 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/pai-rbp1-serbp1-serpine1-monoclonal-mouse-antibody-serbp1-3496/>