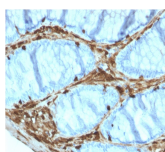


# Annexin A1 Monoclonal Mouse Antibody (ANXA1/3566)



## Product Description

Annexin A1 belongs to the annexin family. It plays a role in glucocorticoid-mediated down-regulation of the early phase of the inflammatory response. In resting conditions Annexin A1 is localized to the cytoplasm. Upon activation it is mobilized to the membrane and secreted. Annexin A1 and contains 4 annexin repeats. A pair of annexin repeats may form one binding site for calcium and a phospholipid. Expression of the ANXA1 gene is upregulated in hairy cell leukemia (HCL). Detection of ANXA1 provides a simple, highly sensitive and specific assay for diagnosing HCL. Annexin A1 has also been found to be protective against DNA damage induced by heat in breast cancer cells, suggesting it is involved in tumor suppressive and protective activities, and also is associated with treatment resistance.

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.

**Catalog number key for antibody number 3566, Annexin A1 / Monoclonal Mouse Antibody (ANXA1/3566)**

## Product attributes

research-areas	Immunology
Antibody number	#3566
Antibody reactivity (target)	Annexin A1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ANXA1/3566
Isotype	IgG2a, kappa
Molecular weight of antigen	38 kDa
Synonyms	Annexin-1; Calpactin-2; Chromobindin-9; Phospholipase A2 inhibitory protein; LPC1; p35
Human gene symbol	ANXA1
Entrez gene ID	301
SwissProt	P04083
Unigene	494173
Immunogen	Recombinant full-length human Annexin A1 protein
Verified antibody applications	Flow (intracellular) (verified), IF (verified), IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Plasma membrane, Nucleus, Secreted (extracellular)
Species reactivity	Human
Positive control	HeLa, A549, A431 and K562 cell lysates, human esophagus, placenta and prostate tissues.
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
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

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>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on June 11, 2026 at 09:53:10 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/annexin-a1-monoclonal-mouse-antibody-anxa1-3566/>