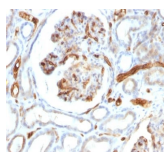


Adiponectin Monoclonal Mouse Antibody (ADPN/1370)



Product Description

Adiponectin is an adipocytokine, hormones produced in adipose tissue. Adiponectin is abundantly present in plasma and has insulin like effect on glucose levels in the blood. Plasma adiponectin levels are low in insulin resistant patients who are obese, have diabetes mellitus type 2 or HIV-lipodystrophy. In women adiponectin levels tend to be higher than in men, which may be due to androgens suppressing adiponectin levels. Furthermore adiponectin and leptin are both indicated in regulating body weight through direct action on the hypothalamus, influencing appetite. Obese people have low adiponectin levels while levels in anorexia patients are high. Adiponectin acts as ligand for various receptors, two of which have been identified, one probably involved in carbohydrate assimilation, the other in tuning the rate of metabolism.

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.

Catalog number key for antibody number 1370, Anti-Adiponectin (ADPN/1370)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#1370
research-areas	Endocrinology, Metabolism
Antibody reactivity (target)	Adiponectin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ADPN/1370
Isotype	IgG2b
Molecular weight of antigen	26-30 kDa
Synonyms	ADPN; ACRP30; Adipocyte Complement-Related Protein; Gelatin binding protein
Human gene symbol	ADIPOQ
Entrez gene ID	9370 (Human)
SwissProt	Q15848 (Human)
Unigene	80485 (Human)
Immunogen	Recombinant human Adiponectin protein fragment
Antibody target cellular localization	Secreted (extracellular)
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL. Immunohistology (formalin) 1-2 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, ELISA For coating, order Ab without BSA, Optimal dilution for a specific application should be determined by user
Positive control	HeLa cells or Kidney
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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
Antibody research areas	Endocrinology, Metabolism
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	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
BNC74	CF®740	742/767	633-685	775/50	CF®740 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 Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on June 7, 2026 at 02:21:38 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mouse-monoclonal-anti-adiponectin-adpn1370/>