

# **TRAcP Monoclonal Mouse Antibody** (ACP5/1070)



## **Product Description**

This antibody recognizes a protein of 35 kDa, which is identified as tartrate-resistant acid phosphatase (TRACP). It exists as two isoforms (5a and 5b). This MAb reacts with both the isoforms. Serum TRACP 5a is secreted by macrophages and dendritic cells and increased in many patients of rheumatoid arthritis.Serum TRAcP 5b is produced from osteoclasts and elevated during bone resorption. TRAcP is an iron containing glycoprotein, which catalyzes the conversion of orthophosphoric monoester to alcohol and orthophosphate. It is the most basic of the acid phosphatases and is the only form not inhibited by L( )-tartrate. TRACP is synthesized as a latent proenzyme and is activated by proteolytic cleavage and reduction. Normally, TRAcP is highly expressed by osteoclasts, activated macrophages, neurons and endometrium during pregnancy. Expression of TRAcP is increased in certain pathological conditions such as Leukemic Reticuloendotheliosis (Hairy Cell Leukemia), Gaucher's Disease, HIV-induced Encephalopathy, Osteoclastoma and in osteoporosis and metabolic bone diseases. Anti-TRACP antibody labels the cells of Hairy Cell Leukemia (HCL) with a high degree of sensitivity and specificity. Other cells stained with this antibody are tissue macrophages and osteoclasts.

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 1070, Anti-TRAcP (ACP5/1070)

#### Call us : 800-304-5357

#### Email: techsupport@biotium.com

### Product attributes

Antibody number	#1070			
Antibody reactivity (target)	TRAcP			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	ACP5/1070			
Isotype	IgG2b, kappa			
Molecular weight	35 kDa			
Synonyms	Acid phosphatase 5, tartrate resistant, ACP5, serum band 5 tartrate-resistant acid phosphatase, SPENCDI, Tartrate-resistant acid ATPase, Tartrate-resistant acid phosphatase type 5, TR-AP, TRACP, TRACP5, TRAP, TrATPase, Type 5 acid phosphatase			
Human gene symbol	ACP5			
Entrez gene ID	54			
SwissProt	P13686			
Unigene	1211			
Immunogen	Recombinant full-length human ACP5 protein			
Verified antibody applications	IHC (FFPE) (verified)			
Antibody target cellular localization	Lysosomes			
Species reactivity	Human, Mouse, Rat			
Species reactivity Antibody application notes	Human, Mouse, Rat Higher concentration may be required for direct detection using primary antibody, roonjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 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 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user			
Species reactivity Antibody application notes Positive control	Human, Mouse, Rat Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 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 minutes, Flow Cytometry 0.5-1 ug/mllion cells/0.1 mL, Optimal dilution for a specific application should be determined by user HepG2, 293T, K562 or RPMI-8226 Cells. Spleen from Hairy Cell Leukemia (HCL) patient.			
Species reactivity Antibody application notes Positive control Shipping condition	Human, Mouse, Rat Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires bolling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/mllion cells/0.1 mL, Optimal dilution for a specific application should be determined by user HepG2, 293T, K562 or RPMI-8226 Cells. Spleen from Hairy Cell Leukemia (HCL) patient. Room temperature			
Species reactivity Antibody application notes Positive control Shipping condition Storage Conditions	Human, Mouse, Rat Higher concentration may be required for direct detection using primary antibody, roonjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 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 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user HepG2, 293T, K562 or RPMI-8226 Cells. Spleen from Hairy Cell Leukemia (HCL) patient. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C			
Species reactivity Antibody application notes Positive control Shipping condition Storage Conditions Shelf life	Human, Mouse, Rat   Higher concentration may be required for direct detection using primary antibody, conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 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 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user   HepG2, 293T, K562 or RPMI-8226 Cells. Spleen from Hairy Cell Leukemia (HCL) patient.   Room temperature   Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C   Guaranteed for at least 24 months from date of receipt when stored as recommended			
Species reactivity Antibody application notes Antibody application notes Antibody application notes Positive control Shipping condition Storage Conditions Shelf life Regulatory status	Human, Mouse, Rat   Higher concentration may be required for direct detection using primary antibody, conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 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 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user   HepG2, 293T, K562 or RPMI-8226 Cells. Spleen from Hairy Cell Leukemia (HCL) patient.   Room temperature   Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C   Guaranteed for at least 24 months from date of receipt when stored as recommended   For research use only (RUO)			
Species reactivity Antibody application notes Antibody application notes Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Human, Mouse, Rat   Higher concentration may be required for direct detection using primary antibody, roonjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 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 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user   HepG2, 293T, K562 or RPMI-8226 Cells. Spleen from Hairy Cell Leukemia (HCL) patient.   Room temperature   Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C   Guaranteed for at least 24 months from date of receipt when stored as recommended   For research use only (RUO)   Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.5% azide, Purified: 0.2 mg/mL in PBS/0.05% BSA, 0.5%			
Species reactivity Antibody application notes Antibody application notes Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	Human, Mouse, Rat   Higher concentration may be required for direct detection using primary antibody, roonjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 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 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user   HepG2, 293T, K562 or RPMI-8226 Cells. Spleen from Hairy Cell Leukemia (HCL) patient.   Room temperature   Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C   Guaranteed for at least 24 months from date of receipt when stored as recommended   For research use only (RUO)   Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.5% azide, Purified: 0.2 mg/mL in PBS/0.5% BSA, 0.5% BSA, prime in PBS/0.5% BSA, 0.5%			
Species reactivity Antibody application notes Antibody application notes Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation Antibody research areas Product origin	Human, Mouse, Rat   Higher concentration may be required for direct detection using primary antibody, roonjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Immunohistology formalin-fixed 0.5-1 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 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution or a specific application should be determined by user   HepG2, 293T, K562 or RPMI-8226 Cells. Spleen from Hairy Cell Leukemia (HCL) patient.   Room temperature   Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C   Guaranteed for at least 24 months from date of receipt when stored as recommended   For research use only (RUO)   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, Purified: 0.2 mg/mL in PBS/0.05% azide, PMP conjugates: 0.1 mg/mL in PBS/0.05% azide, PMP time PBS without azide   Metabolism   Product may contain either bovine serum albumin (BSA) from bovine serum (BSA tures), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.			

Cell/tissue expression

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 July 13, 2025 at 12:21:33 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-tracp-tartarate-resistant-acid-phosphatase-acp51070/