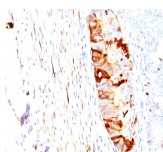


Aldehyde Dehydrogenase 1 Monoclonal Mouse Antibody (ALDH1A1/1382)



Product Description

Aldehyde dehydrogenase 1 family member A1 (ALDH1A1), also known as retinal dehydrogenase 1, belongs to the aldehyde dehydrogenase enzyme family, which is involved in the metabolism of alcohol. ALDH1A1 is predominantly expressed in the epithelium of testis, brain, eye, liver, kidney, as well as neural and hematopoietic stem cells. Reportedly, high ALDH1A1 expression is found in solitary fibrous tumor (SFT) and hemangiopericytoma (HPC), compared to meningiomas and synovial sarcomas. In combination with CD34, ALDH1A1 may be useful for the differentiation among SFT, HPC, meningioma, and synovial sarcoma.

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 1382, ALDH1A1 Monoclonal Mouse Antibody (ALDH1A1/1382)

Product attributes

research-areas	Metabolism
Antibody number	#1382
Antibody reactivity (target)	Aldehyde dehydrogenase 1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	ALDH1A1/1382
Isotype	IgG
Molecular weight	56 kDa
Synonyms	Aldehyde dehydrogenase family 1 member A1; ALDH1A1; Retinal dehydrogenase 1; Retinaldehyde Dehydrogenase 1; Epididymis Luminal Protein
Human gene symbol	ALDH1A1
Entrez gene ID	216
SwissProt	P00352
Unigene	76392
Immunogen	Purified recombinant fragment of human ALDH1A1 (aa 315-434), expressed in HK293 Cells.
Antibody target cellular localization	Cytoplasmic
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	HepG2 or A549 cells. Colon or Testicular Carcinoma.
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
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.

References

Xu J, Muller S, et al. Eur J Cancer. 2012 Jul;48(11):1682-91

This datasheet was generated on August 23, 2025 at 09:56:04 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/aldh1a1-monoclonal-mouse-antibody-aldh1a1-1382/>