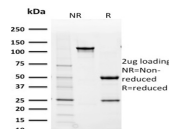


Aldo-keto Reductase Family 1 Member B1 Monoclonal Mouse Antibody (CPTC-AKR1B1-2)



Product Description

Aldo-keto Reductase Family 1 Member B1 (AKR1B1), also known as aldose reductase or aldehyde reductase, is a member of the aldo/keto reductase superfamily, which consists of more than 40 known enzymes and proteins. This protein catalyzes the reduction of a number of aldehydes, including the aldehyde form of glucose, and is thereby implicated in the development of diabetic complications by catalyzing the reduction of glucose to sorbitol. It has also been shown to have decreased expression in adrenocortical cancer, and possibly play a role in adrenal tumorigenesis. It has been suggested that AKR1B1 could be investigated as a marker of malignancy for adrenal tumor diagnosis. 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 2592, Aldo-keto Reductase Family 1 Member B1 Monoclonal Mouse Antibody (CPTC-AKR1B1-2)**

Product attributes

research-areas	Metabolism
Antibody number	#2592
Antibody reactivity (target)	Aldo-keto Reductase Family 1 Member B1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-AKR1B1-2
Isotype	IgG1, kappa
Molecular weight	37 kDa
Synonyms	Aldo-keto reductase family 1 member B1; AKR1B1; Aldehyde Reductase 1; Aldose reductase; Lii5-2 CTCL tumor antigen; low Km aldose reductase
Human gene symbol	AKR1B1
Entrez gene ID	231
SwissProt	P15121
Unigene	521212
Immunogen	Recombinant full-length human AKR1B1 protein
Verified antibody applications	IF (verified), IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Cytoplasmic
Species reactivity	Human
Positive control	549, 293T, A431, HeLa, HepG2, MOLT4 and Raji whole cell lysates; HeLa cells; Human colon carcinoma tissue.
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
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

Dye Features

[CF®405S Features](#)
[CF®488A Features](#)
[CF®568 Features](#)
[CF®594 Features](#)
[CF®640R Features](#)
[CF®647 Features](#)

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 May 18, 2024 at 05:16:59 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/aldo-keto-reductase-family-1-member-b1-monoclonal-mouse-antibody-cptc-akr1b1-2/>