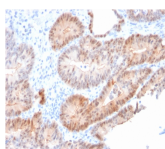


Aldo-keto Reductase Family 1 Member C2 Monoclonal Mouse Antibody (CPTC-AKR1C2-1)



Product Description

Aldo-keto Reductase Family 1 Member C2 (AKR1C2), also known as Dihydrodiol Dehydrogenase 2 (DDH2), is a member of the aldo/keto reductase superfamily, which consists of more than 40 known enzymes and proteins. These enzymes catalyze the conversion of aldehydes and ketones to their corresponding alcohols by utilizing NADH and/or NADPH as cofactors. The enzymes display overlapping but distinct substrate specificity. This enzyme catalyzes the reaction of progesterone to the inactive form 20-alpha-hydroxy-progesterone.

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 2255, Aldo-keto Reductase Family 1 Member C2 / DD2 Monoclonal Mouse Antibody (CPTC-AKR1C2-1)

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

Product attributes

research-areas	Metabolism
Antibody number	#2255
Antibody reactivity (target)	Aldo-keto Reductase Family 1 Member C2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-AKR1C2-1
Isotype	IgG2a, kappa
Molecular weight	37 kDa
Synonyms	Aldo-keto Reductase Family 1 Member C2; AKR1C2; 3-alpha-HSD3; AKR1C-pseudo protein; BAP; Dihydrodiol dehydrogenase 2; DD-2; MCDR2; HAKRD; DD; DDH2; HBAB; Pseudo-chlordecone reductase; SRXY8
Human gene symbol	AKR1C2
Entrez gene ID	1646
SwissProt	P52895
Unigene	567256
Immunogen	Recombinant human full-length AKR1C2 protein
Verified antibody applications	IHC (FFPE), WB (verified)
Antibody target cellular localization	Cytoplasmic
Species reactivity	Human
Positive control	HeLa, K-562, A431, HepG2, A549 cells. Human liver or stomach 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
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 January 5, 2026 at 07:10:32 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/aldo-keto-reductase-family-1-member-c2-dd2-monoclonal-mouse-antibody-cptc-akr1c2-1/>