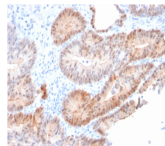


Aldo-keto Reductase Family 1 Member C2 Monoclonal Mouse Antibody (CPTC-AKR1C2-1)



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)**

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; BABP; Dihydrodiol dehydrogenase 2; DD-2; MCDR2; HAKRD; DD; DDH2; HBAB; Pseudo-chlorocone 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) (verified), 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

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
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, LI-COR Bioscience.