## **UGT1A9 Monoclonal Mouse Antibody** (**UGT1A9**/1229)

## **Product Description**

This antibody recognizes a protein of about 60 kDa, which is identified as human UGT1A9. It does not cross-react with the other UGT1A isoforms including UGT1A7, UGT1A8, and UGT1A10 and shows a high degree of specificity. UGT1A9 is a UDP-glucuronosyltransferase, an enzyme of the glucuronidation pathway that transforms small lipophilic molecules, such as steroids, bilirubin, hormones, and drugs, into water-soluble, excretable metabolites. This MAb binds to human hepatocytes and the majority of human hepatocellular carcinomas (HCC s). In frozen sections, it stains hepatic cells and may be used as a marker of the liver. This MAb also binds to cell preparations of hepatocellular carcinoma biopsies or cell lines

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 <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1229, Anti-UGT1A9 (UGT1A9/1229)

## Product attributes

Product attributes				
Antibody number	#1229			
Antibody reactivity (target)	UGT1A9			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	UGT1A9/1229			
Isotype	IgG2b, kappa			
Molecular weight	60 kDa			
Synonyms	GNT1; HLUGP4; lugP4; UDP-glucuronosyltransferase 1-l; UDP-glucuronosyltransferase 1A9 (UGT1A9); UDPGT 1-9; UDPGT; UGT-1I; UGT1*9; UGT1-09; UGT1.9; Ugt1a9; UGT1I			
Human gene symbol	UGT1A9			
Entrez gene ID	54600			
SwissProt	O60656			
Unigene	554822			
Immunogen	A synthetic peptide specific to human UGT1A9 protein			
Antibody target cellular localization	Endoplasmic reticulum, Microsomes			
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 cells. Liver or Hepatocellular Carcinoma (HCC)			
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			
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended			
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% BSA-free: 1 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide			
	Metabolism			
Antibody research areas	Wetabolish			
Antibody research areas 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.			

Call us: 800-304-5357 Email: btinfo@biotium.com

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

Shingo Oda, Miki Nakajima, Masahiko Hatakeyama, Tatsuki Fukami, and Tsuyoshi Yokoi. Preparation of a Specific Monoclonal Antibody against Human UDP-Glucuronosyltransferase (UGT) 1A9 and Evaluation of UGT1A9 Protein Levels in Human Tissues. Drug Metabolism and Disposition, 40:1620-1627, 2012

This datasheet was generated on November 5, 2025 at 08:18:55 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/monoclonal-anti-ugt1a9-udp-glucuronosyltransferase-1-family-polypeptide-a9-ugt1a9133/">https://biotium.com/product/monoclonal-anti-ugt1a9-udp-glucuronosyltransferase-1-family-polypeptide-a9-ugt1a9133/</a>