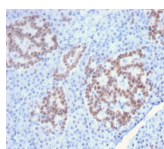


HNF1A Monoclonal Mouse Antibody (HNF1A/2087)



Product Description

HNF1A belongs to the homeobox protein family and is an essential transcription factor for many hepatic genes involved in detoxification, homeostasis and metabolisms of glucose, lipid, steroid and amino acid. In addition, HNF1A is an important component of the transcriptional networks governing embryonic pancreas development and differentiation. as well as maintaining the growth and function of islet cells in adult. HNF1A (Hepatocyte nuclear factor 1 alpha) is a transcription factor that is known to regulate pancreatic differentiation and maintain homeostasis of endocrine pancreas. Recently, genome-wide association studies have implicated HNF1A as a susceptibility gene for pancreatic cancer.

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 2087, Anti-HNF1A (HNF1A/2087)

Product attributes

Antibody number	#2087
Antibody reactivity (target)	HNF1A
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	HNF1A/2087
Isotype	IgG1, kappa
Molecular weight	79 kDa
Synonyms	Albumin proximal factor; Hepatic nuclear factor 1 alpha; Hepatic transcription factor 1 alpha; HNF-1A; Interferon production regulator factor; LFB1; LFB1 hepatic nuclear factor; Liver specific transcription factor LFB1; Maturity onset diabetes of the young 3 (MODY3); TCF1; Transcription factor 1 hepatic
Human gene symbol	HNF1A
Entrez gene ID	6927
SwissProt	P20823
Unigene	654455
Immunogen	Recombinant fragment (around aa 214-339) of human HNF1A protein (exact sequence is proprietary)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Positive control	Jurkat cells. Small Intestine or Pancreas.
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 April 18, 2025 at 11:40:24 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/hnf1a-monoclonal-mouse-antibody-hnf1a-2087/>