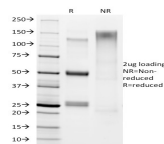


Helicobacter pylori Monoclonal Mouse Antibody (HP/1336)



Product Description

The spiral shaped bacterium *Helicobacter pylori* is strongly associated with inflammation of the stomach and is also implicated in the development of gastric malignancy. *H. pylori* is known to cause peptic ulcers and chronic gastritis in human. It is associated with duodenal ulcers and may be involved in development of adenocarcinoma and low-grade lymphoma of mucosa associated lymphoid tissue in the stomach. This antibody stains the individual *H. pylori* bacterium when it presents on the surface of the epithelium or in the cytoplasm of the epithelial cells in biopsy tissue sections from the antrum and body of the stomach.

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 1336, Anti-Helicobacter pylori (HP/1336)

Product attributes

Antibody number	#1336
Antibody reactivity (target)	<i>Helicobacter pylori</i>
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	HP/1336
Isotype	IgG1, kappa
Molecular weight of antigen	Not Known
Synonyms	<i>H. pylori</i>
Entrez gene ID	Not Known
SwissProt	Not Known
Unigene	Not Known
Immunogen	Total sonicate of <i>Helicobacter pylori</i>
Species reactivity	<i>Helicobacter pylori</i>
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	<i>Helicobacter pylori</i> infected stomach biopsy
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
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 (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Antibody research areas	Infectious disease, Microbiology

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 June 15, 2026 at 02:03:03 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/helicobacter-pylori-monoclonal-mouse-antibody-hp-1336/>