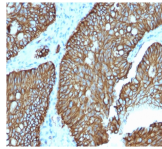


Cadherin 17 / LI Cadherin Monoclonal Mouse Antibody (CDH17/2618)



Product Description

Cadherin 17, also known as LI cadherin (for liver-intestine cadherin), is a member of the cadherin family of calcium-dependent adhesion molecules. Cadherins mediate cell-cell interactions via hemophilic binding, and are critical to the maintenance of tissue structure and morphogenesis. Cadherins each contain a large extracellular domain at the amino terminus, which is characterized by a series of five homologous repeats, the most distal of which is thought to be responsible for binding specificity. The relatively short carboxy terminal, intracellular domain interacts with a variety of cytoplasmic proteins, including beta-catenin, to regulate cadherin function. Cadherin 17 expression is restricted to liver and intestine tissues and is specifically localized to the basolateral domain of hepatocytes and enterocytes.

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 2618, Cadherin 17 / LI Cadherin Monoclonal Mouse Antibody (CDH17/2618)

Product attributes

Antibody number	#2618
Antibody reactivity (target)	Cadherin 17, LI Cadherin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CDH17/2618
Isotype	IgG2b, kappa
Molecular weight	120 kDa
Synonyms	Cadherin 17; Cadherin-17; CDH17; LI-cadherin; Liver-intestine cadherin; human peptide transporter 1 (HPT-1); Intestinal peptide-associated transporter HPT-1
Human gene symbol	CDH17
Entrez gene ID	1015
SwissProt	Q12864
Unigene	591853
Immunogen	Recombinant fragment (around aa 242-418) of human Cadherin 17 protein (CDH17) (exact sequence is proprietary)
Verified antibody applications	Flow (verified), IHC (FFPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Positive control	MCF-7, HepG2 or HT29 cells. Stomach, Small Intestine, Colon or liver.
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 research areas	Cell adhesion
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.
Cell/tissue expression	Intestine, Liver

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 July 13, 2025 at 12:13:10 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cadherin-17-li-cadherin-monoclonal-mouse-antibody-cdh17-2618/>