

Galectin-13 / Placental Protein 13 Monoclonal Mouse Antibody (PP13/1162)

Product Description

This antibody recognizes a 32 kDa protein, which is identified as homodimer of galectin-13 (also known as PP13). Galectins are a family of soluble beta-galactoside-binding lectins that modulate cell-to-cell adhesion and cell-to-extracellular matrix (ECM) interactions and play a role in tumor progression, pre-mRNA splicing and apoptosis. One such member, galectin-13, is a 139 amino acid protein that contains one galectin domain. Possessing lysophospholipase activity, galectin-13 exists as a disulfide-linked homodimer. Galectin-13 is suggested to have a developmental role in the placenta and may display immunobiological activity in feto-maternal blood-spaces. Screening for galectin-13 in maternal serum during the first trimester of pregnancy may serve as a diagnostic marker to predict preterm pre-eclampsia.

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 1162, Anti-Galectin-13 (PP13/1162)

Product attributes

Antibody number	#1162
Antibody reactivity (target)	Galectin-13, Placental Protein 13
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	PP13/1162
Isotype	IgG2b, kappa
Molecular weight of antigen	16 kDa (monomer); 32 kDa (homodimer)
Synonyms	Gal-13; GAL13; Galactoside-binding soluble lectin 13; Galectin-13; Galectin13; Lectin galactoside binding soluble 13; LGALS13; PLAC8; Placental protein 13 (PP13); Placental tissue protein 13
Human gene symbol	LGALS13
Entrez gene ID	29124
SwissProt	Q9UHV8
Unigene	23671
Immunogen	Recombinant human Galectin-13 protein fragment (aa23-134) (exact sequence is proprietary)
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 0.5-1 ug/mL. Immunohistology (formalin): 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris, 1 mM EDTA pH 9.0 or 10 mM citrate buffer pH 6.0 for 10-20 min followed by cooling at RT for 20 min. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. ELISA For coating, order Ab without BSA. Optimal dilution for a specific application should be determined by user
Positive control	K-562 cells. Placenta or Spleen.
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% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Antibody research areas	Developmental biology, Immunology
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 # 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 13, 2026 at 08:43:54 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mouse-monoclonal-anti-galectin-13-placental-protein-13-pp131162/>