## Galectin-13 / Placental Protein 13 Monoclonal Mouse Antibody (PP13/1161)

## **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 <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 1161, Anti-Galectin-13 (PP13/1161)

## Product attributes

Product attributes				
Antibody number	#1161			
Antibody reactivity (target)	Galectin-13, Placental Protein 13			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	PP13/1161			
Isotype	IgG2b, kappa			
Molecular weight	16 kDa (monomer); 32 kDa (homodimer)			
Synonyms	Gal-13; GAL13; Galactoside-binding soluble lectin 13; Galectin 13; Galectin 13; Lectin galactoside binding soluble 13; LGALS13; PLAC8; Placental protein 13 (PP13); Placental tissu 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 (FFPE) (verified), WB (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, Western blotting 0.5-1 ug/mL, 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/m in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL PBS without azide			
Antibody research areas	Developmental biology, Immunology			
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: techsupport@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.

This datasheet was generated on October 27, 2025 at 08:58:51 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/mouse-monoclonal-anti-galectin-13-placental-protein-13-pp131161/">https://biotium.com/product/mouse-monoclonal-anti-galectin-13-placental-protein-13-pp131161/</a>