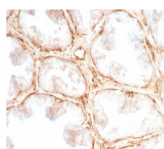


Galectin-1 / Human Placental Lactogen Recombinant Monoclonal Rabbit Antibody (GAL1/2499R)



Product Description

Galectin-1 is a member of the beta-galactoside-binding family and is a dimeric protein of 14 kD participating in a variety of normal and pathological processes, including cancer progression. Galectin-1 can affect the proliferation of normal and malignant cells. Inhibition of cell growth is observed in a lactose-dependent manner as lower concentrations of the lectin stimulate cell proliferation. Galectin-1 may also be implicated in the induction of apoptosis of activated T cells through the binding of exogenous galectin-1 to CD45 molecules present on the surface of lymphocytes. Galectin-1, reported to be present either at the surface of cancer cells or accumulated around these cells could act as an immunological shield to protect against a T cell immune response and provide an advantage for survival.

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 2499, Anti-Galectin-1 / Human Placental Lactogen (GAL1/2499R)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#2499
Antibody reactivity (target)	Galectin-1, Human Placental Lactogen
Antibody type	Primary
Host species	Rabbit
Clonality	Monoclonal, Recombinant Monoclonal
Clone	GAL1/2499R
Isotype	IgG, kappa
Molecular weight of antigen	14 kDa
Synonyms	14kDa laminin-binding protein; Beta-galactoside-binding lectin L-14-I; Gal-1; GAL1; Galaptin; Galectin-1; HLBP14; HPL; Lactose-binding lectin 1; Lect14; Lectin galactoside-binding soluble 1; LGALS1; MAPK activating protein MP12; S-Lac lectin 1
Human gene symbol	LGALS1
Entrez gene ID	3956
SwissProt	P09382
Unigene	445351
Immunogen	Recombinant human Galectin-1 protein fragment (aa12-108) (exact sequence is proprietary)
Verified antibody applications	IHC (FPPE) (verified), WB (verified)
Antibody target cellular localization	Extracellular matrix
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, ELISA: 2-4 ug/m for coating order Ab without BSA, Immunohistology (formalin): 0.5-1 ug/mL for 30 minutes at RT, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user
Positive control	HeLa, K562 or 293 cells. Prostate, Kidney, Placenta, Stomach, Skin, Spleen, Brain or Heart.
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	Cancer, 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 15, 2026 at 02:05:29 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/galectin-1-human-placental-lactogen-recombinant-monoclonal-rabbit-antibody-gal1-2499r/>