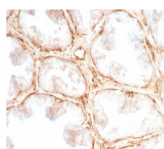


# Galectin-1 / Human Placental Lactogen Recombinant Monoclonal Rabbit Antibody (GAL1/2499R)



## 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](mailto: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

<b>Antibody number</b>	#2499
<b>Antibody reactivity (target)</b>	Galectin-1, Human Placental Lactogen
<b>Antibody type</b>	Primary
<b>Host species</b>	Rabbit
<b>Clonality</b>	Monoclonal, Recombinant Monoclonal
<b>Clone</b>	GAL1/2499R
<b>Isotype</b>	IgG
<b>Molecular weight of antigen</b>	14 kDa
<b>Synonyms</b>	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
<b>Human gene symbol</b>	LGALS1
<b>Entrez gene ID</b>	3956
<b>SwissProt</b>	P09382
<b>Unigene</b>	445351
<b>Immunogen</b>	Recombinant human Galectin-1 protein fragment (aa12-108) (exact sequence is proprietary)
<b>Verified antibody applications</b>	IHC (FFPE) (verified), WB (verified)
<b>Antibody target cellular localization</b>	Extracellular matrix
<b>Species reactivity</b>	Human
<b>Antibody application notes</b>	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
<b>Positive control</b>	HeLa, K562 or 293 cells. Prostate, Kidney, Placenta, Stomach, Skin, Spleen, Brain or Heart.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Cancer, Immunology
<b>Product origin</b>	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	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 Features</a>
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 8, 2026 at 09:24:37 PM. 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/>