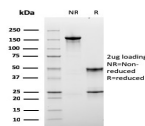


Lymphocyte Activation Gene 3 Monoclonal Mouse Antibody (LAG3/3261)



Product Description

LAG-3 (also called CD223) is a high affinity MHC class II ligand present on the surface of CD4 CD8 T cells and NK cell, with shared homology in structure to CD4 molecules. It has a glutamic acid-proline (EP) repetitive sequence found in other functionally distinct mammalian, parasitic, and bacterial proteins that may influence a conserved biological function. LAG-3 CD4 CD8 T cells can associate with the T cell receptor (TCR) and downregulate TCR signaling in vitro. LAG-3 inhibits CD4-dependent T cell function via its cytoplasmic domain. LAG-3 Lys-468 within a conserved 'KIEELE' motif is essential for interaction with downstream signaling molecules. Furthermore, as a checkpoint inhibitor target, it may be superior to CTLA-4 and PD-1 since both antibodies only activate effector T-cells, whereas an antagonist LAG-3 antibody can both activate T effector cells (by downregulating the LAG-3 inhibiting signal into pre-activated LAG-3 cells) and inhibit induced (i. e. antigen-specific) Treg suppressive activity.

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 3261, Anti-Lymphocyte Activation Gene 3 (LAG3/3261)

Product attributes

Antibody number	#3261
Antibody reactivity (target)	Lymphocyte Activation Gene 3
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	LAG3/3261
Isotype	IgG2a, kappa
Molecular weight of antigen	70 kDa
Synonyms	CD223; FDC protein; LAG-3
Human gene symbol	LAG3
Entrez gene ID	3902
SwissProt	P18627
Unigene	409523
Immunogen	Recombinant full-length human LAG3 protein
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Human tonsil or Hodgkin's lymphoma (IHC-P).
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
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	NK cells, T-cells

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 21, 2026 at 04:18:55 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/lymphocyte-activation-gene-3-monoclonal-mouse-antibody-lag3-3261/>