

PRODUCT INFORMATION



Galectin 9 Short Variant (human, recombinant)

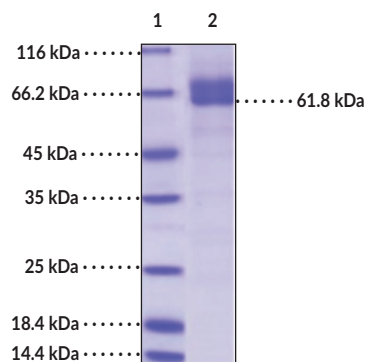
Item No. 32012

Overview and Properties

Synonyms: Gal-9S, HUAT, Human Urate Transporter, Lectin Galactoside-binding Soluble 9, LGALS9A
Source: Recombinant N-terminal human IgG1 Fc-tagged galectin 9 expressed in HEK293 cells
Amino Acids: 2-323 (full length)
Molecular Weight: 61.8 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥90% estimated by SDS-PAGE
Supplied in: Lyophilized from sterile PBS, pH 7.4
Endotoxin Testing: < 1.0 EU/μg, determined by the LAL endotoxin assay

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers

Lane 2: Galectin 9 Short Variant

SDS-PAGE Analysis of Galectin 9 Short Variant.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/18/2020

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Galectin 9 is a β -galactoside-binding protein encoded by the *LGALS9* gene and a member of the galectin family.¹ It contains two non-identical carbohydrate recognition domains (CRDs) joined by a peptide linker, the length of which depends on alternative splicing of the *LGALS9* gene, which produces a short, medium, or long variant of the protein.^{1,2} The non-identical CRDs are responsible for binding to different types of saccharide ligands with the N-terminal CRD (NCRD) preferring more complex glycoconjugates.¹ Galectin 9 is localized to the cytosol, nucleus, and extracellular matrix and is expressed primarily in immune-related tissues and cells.³ When expressed on immune or tumor cells, galectin 9 binds to glycosylated sites on the immunoregulatory protein TIM-3 and activates signaling that impairs immune synapse formation leading to T cell anergy or apoptosis.⁴ It also binds to the transmembrane glycoprotein CD44 on osteoblasts, leading to their proliferation, and it has a role in cell adhesion by preventing hyaluronic acid from binding to CD44.⁵ The short variant of galectin 9 acts as an eosinophil chemoattractant.⁶ It also activates CD4⁺ T cell populations and induces naïve T cells to acquire a central memory T cell (T_{CM}) phenotype.⁷ Galectin 9 was originally identified as a tumor antigen in patients with Hodgkin's disease.¹ It is expressed to a lower extent in tumor cells compared with non-tumor cells but is increased in leukemia and colon cancer cell lines.³ Increased expression of the short variant of galectin 9 inhibits or enhances adhesion in breast or colon cancer cell lines, respectively. The expression of the short variant of galectin 9 is increased by interferon γ (IFN- γ) in synovial fibroblasts isolated from patients with rheumatoid arthritis.⁸ Cayman's Galectin 9 Short Variant (human, recombinant) protein is a disulfide-linked homodimer. The reduced monomer, comprised of galectin 9 (amino acids 2-323) fused to human IgG1 Fc at its N-terminus, consists of 554 amino acids and has a calculated molecular weight of 61.8 kDa.

References

1. Nagae, M., Nishi, N., Murata, T., *et al.* Crystal structure of the galectin-9 N-terminal carbohydrate recognition domain from *Mus musculus* reveals the basic mechanism of carbohydrate recognition. *J. Biol. Chem.* **281(47)**, 35884-35893 (2006).
2. Chabot, S., Kashio, Y., Seki, M., *et al.* Regulation of galectin-9 expression and release in Jurkat T cell line cells. *Glycobiology* **12(2)**, 111-118 (2002).
3. Heusschen, R., Griffioen, A.W., and Thijssen, V.L. Galectin-9 in tumor biology: A jack of multiple trades. *Biochim. Biophys. Acta* **1836(1)**, 177-185 (2013).
4. Wolf, Y., Anderson, A.C., and Kuchroo, V.K. TIM3 comes of age as an inhibitory receptor. *Nat. Rev. Immunol.* **20(3)**, 173-185 (2020).
5. John, S. and Mishra, R. Galectin-9: From cell biology to complex disease dynamics. *J. Biosci.* **41(3)**, 507-534 (2016).
6. Sato, M., Nishi, N., Shoji, H., *et al.* Functional analysis of the carbohydrate recognition domains and a linker peptide of galectin-9 as to eosinophil chemoattractant activity. *Glycobiology* **12(3)**, 191-197 (2002).
7. Gooden, M.J.M., Wiersma, V.R., Samplonius, D.F., *et al.* Galectin-9 activates and expands human T-helper 1 cells. *PLoS One* **8(5)**, e65616 (2013).
8. Pearson, M.J., Bik, M.A., Ospelt, C., *et al.* Endogenous galectin-9 suppresses apoptosis in human rheumatoid arthritis synovial fibroblasts. *Sci. Rep.* **8(1)**, 12887 (2018).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM