

PRODUCT INFORMATION



VISTA (human, recombinant)

Item No. 32086

Overview and Properties

Synonyms: DD1 α , Death Domain 1 α , PD-1H, Programmed Cell Death 1 Homolog, SISP1, Stress-Induced Secreted Protein 1, V-domain Ig Suppressor of T cell activation, V-set Domain-Containing Immunoregulatory Receptor, V-set Immunoregulatory Receptor

Source: Recombinant human C-terminal His-tagged VISTA expressed in HEK293 cells

Amino Acids: 33-194

Uniprot No.: Q9H7M9

Molecular Weight: 19.6 kDa

Storage: -80°C (as supplied)

Stability: ≥ 1 year

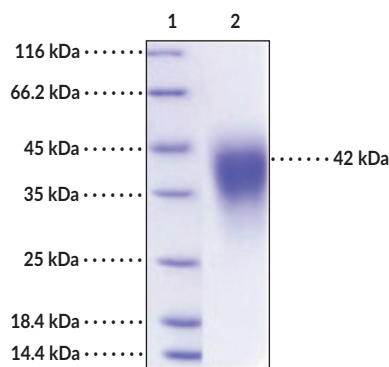
Purity: $\geq 95\%$ estimated by SDS-PAGE

Supplied in: Lyophilized from sterile PBS, pH 7.4

Endotoxin Testing: < 1.0 EU/ μ g, 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: VISTA

SDS-PAGE Analysis of VISTA. This protein has a calculated molecular weight of 19.6 kDa. It has an apparent molecular weight of approximately 38 to 42 kDa by SDS-PAGE under reducing conditions due to apparent post-translational modifications.

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, 06/26/2020

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

V-Domain immunoglobulin (Ig) suppressor of T cell activation (VISTA) is a type I transmembrane receptor and negative checkpoint regulator encoded by *VSIR* in humans.¹ It is composed of an N-terminal extracellular domain, containing a signaling sequence, Ig V domain, and stalk region, a transmembrane domain, and a cytoplasmic tail that contains a conserved Src homology 2 (SH2) binding motif. It is expressed in myeloid cells, monocytes, macrophages, dendritic cells, naïve CD4⁺ and FoxP3⁺ regulatory T cells, and natural killer (NK) cells.² VISTA arrests cell division and suppresses early T cell receptor activation when expressed on CD4⁺ T cells, and suppresses antigen-specific T cell activation when expressed on antigen presenting cells.³ *Vsir*^{-/-} mice are resistant to tumor induction compared with wild-type mice in a GL261 mouse glioma model.^{2,3} Monocyte levels of VISTA are increased in patients with HIV compared with healthy controls.⁴ *VSIR* expression is increased in the circulating immune cells of patients with systemic lupus erythematosus (SLE) and in skin biopsies from patients with discoid lupus erythematosus (DLE). Cayman's VISTA (human, recombinant) protein consists of 173 amino acids, has a calculated molecular mass of 19.6 kDa, and a predicted N-terminus of Phe33 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the molecular mass of the protein is 38-42 kDa due to apparent post-translational modifications.

References

1. Nowak, E.C., Lines, J.L., Varn, F.S., *et al.* Immunoregulatory functions of VISTA. *Immunol. Rev.* **276**(1), 66-79 (2017).
2. ElTanbouly, M.A., Croteau, W., Noelle, R.J., *et al.* VISTA: A novel immunotherapy target for normalizing innate and adaptive immunity. *Semin. Immunol.* **42**, 101308 (2019).
3. Ni, L. and Dong, C. New checkpoints in cancer immunotherapy. *Immunol. Rev.* **276**(1), 52-65 (2017).
4. Wang, G., Tai, R., Wu, Y., *et al.* The expression and immunoregulation of immune checkpoint molecule VISTA in autoimmune diseases and cancers. *Cytokine Growth Factor Rev.* **52**, 1-14 (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