

PRODUCT INFORMATION

JAK3 JH2 Domain (human, recombinant; aa 511-790) - Biotinylated
Item No. 42276

Overview and Properties

Synonyms: Janus-associated Kinase 3, Leukocyte Janus Kinase, L-JAK, Tyrosine-protein Kinase JAK3
Source: Active recombinant human C-terminal AVI- and His-tagged JAK3 JH2 domain expressed in insect cells (sf9)
Amino Acids: 511-790
Uniprot No.: P52333
Molecular Weight: 34 kDa
Storage: -80°C (as supplied)
Stability: ≥6 months
Purity: ≥90% estimated by SDS-PAGE
Supplied in: 40 mM Tris-HCl, pH 8.0, with 110 mM sodium chloride, 2.2 mM potassium chloride, 20% glycerol, and 3 mM DTT

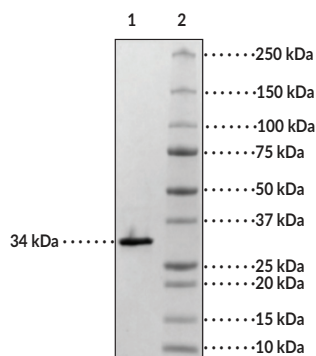
Protein

Concentration: *batch specific* mg/ml

Bioactivity: See figures for details

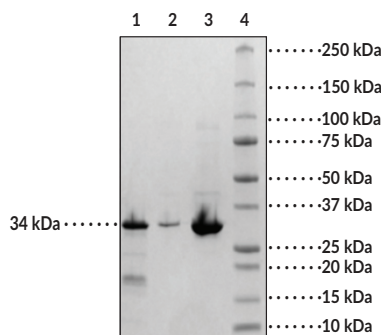
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: JAK3 JH2 Domain
Lane 2: MW Markers

SDS-PAGE Analysis of JAK3 JH2 Domain.



Lane 1: Beads
Lane 2: Flow Through
Lane 3: Positive Control (JAK3 JH2 Domain)
Lane 4: MW Markers

Biotin-Advin pull down SDS-PAGE Analysis of JAK3 JH2 Domain.

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, 10/29/2024

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

JAK3 is a non-receptor tyrosine kinase that has roles in cytokine signaling and immune cell function.^{1,2} It is composed of N-terminal FERM and SH2 domains, an autoinhibitory JH2 pseudokinase domain, and a C-terminal kinase domain.²⁻⁴ JAK3 is constitutively expressed in natural killer (NK) cells and thymocytes and expressed upon cell activation in T cells, B cells, and monocytes.⁵ Following cytokine binding to the IL-2 receptor (IL-2R), IL-4R, IL-7R, IL-9R, IL-15R, or IL-21R, JAK3 binds to the γ_c subunit of the receptor and induces heterodimerization of the receptor subunits and activation of STAT transcription factors.^{1,2,5} Through activation of these receptors, JAK3-mediated signaling is involved in T cell proliferation, differentiation, and survival, B cell differentiation and function, and macrophage activation, among other activities.^{1,2} Loss-of-function mutations in JAK3 are associated with autosomal recessive severe combined immunodeficiency disease (SCID), while gain-of-function mutations are associated with immune dysregulation and blood cancers, including myeloproliferative neoplasms, T cell lymphomas and leukemias, NK lymphoma-leukemia, and acute lymphoblastic leukemia.^{2,6,7} Cayman's JAK3 JH2 Domain (human, recombinant; aa 511-790) – Biotinylated protein can be used for avidin-binding assays and has a calculated molecular weight of 34 kDa.

References

1. Benczik, M., and Gaffen, S.L. The interleukin (IL)-2 family cytokines: Survival and proliferation signaling pathways in T lymphocytes. *Immunol. Invest.* **33**(2), 109-142 (2004).
2. Liongue, C., Ratnayake, T., Basheer, F., et al. Janus kinase 3 (JAK3): A critical conserved node in immunity disrupted in immune cell cancer and immunodeficiency. *Int. J. Mol. Sci.* **25**(5), 2977 (2024).
3. Leonard, W.J., and O'Shea, J.J. Jaks and STATs: Biological implications. *Annu. Rev. Immunol.* **16**, 293-322 (1998).
4. Shanarinen, P., and Silvennoinen, O. The pseudokinase domain is required for suppression of basal activity of Jak2 and Jak3 tyrosine kinases and for cytokine-inducible activation of signal. *J. Biol. Chem.* **277**(49), 47954-47963 (2002).
5. Notarangelo, L.D., and Candotti, F. JAK3-deficient severe combined immunodeficiency. *Radiol. Clin. North Am.* **20**(1), 97-111 (2000).
6. Ott, N., Faletti, L., Heeg, M., et al. JAKs and STATs from a clinical perspective: Loss-of-function mutations, gain-of-function mutations, and their multidimensional consequences. *J. Clin. Immunol.* **43**(6), 1326-1359 (2023).
7. Philips, R.L., Wang, Y., Cheon, H., et al. The JAK-STAT pathway at 30: Much learned, much more to do. *Cell* **185**(21), 3857-3876 (2022).

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