

PRODUCT INFORMATION

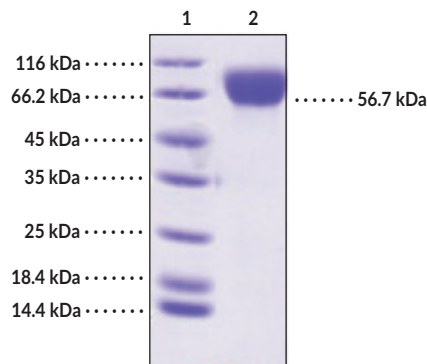


c-Kit/CD117 GNNK- Isoform Extracellular Domain (human, recombinant) Item No. 33980

Overview and Properties

- Synonyms:** c-Kit(-), c-Kit Isoform 2, Cluster of Differentiation 117, Mast/Stem Cell Growth Factor Receptor Kit, Proto-oncogene c-Kit, SCFR, v-Kit Hardy-Zuckerman 4 Feline Sarcoma Viral Oncogene Homolog
- Source:** Active recombinant human C-terminal His-tagged c-Kit GNNK-isoform extracellular domain expressed in HEK293 cells
- Amino Acids:** 26-516
- Uniprot No.:** P10721-2
- Molecular Weight:** 56.7 kDa
- Storage:** -80°C (as supplied)
- Stability:** ≥1 year
- Purity:** ≥97% 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
- 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.*

Image



Lane 1: MW Markers
Lane 2: c-Kit/CD117 GNNK- Isoform Extracellular Domain

SDS-PAGE Analysis of c-Kit/CD117 GNNK- Isoform Extracellular Domain. This protein has a calculated molecular weight of 56.7 kDa. It has an apparent molecular weight of approximately 86 kDa by SDS-PAGE under reducing conditions due to glycosylation.

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, 02/17/2023

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

c-Kit, also known as CD117, is a type III receptor tyrosine kinase involved in the proliferation, differentiation, and survival of hematopoietic cells.^{1,2} It is composed of an N-terminal extracellular domain (ECD) containing five immunoglobulin-like (Ig-like) domains, short transmembrane and juxtamembrane domains, and a C-terminal intracellular region containing the kinase domain. Alternative splicing of *KIT* produces two isoforms, GNNK+ and GNNK-, that contain or do not contain the tetrapeptide GNNK, respectively, in the juxtamembrane region of the ECD.³ c-Kit is expressed in hematopoietic stem and progenitor cells, as well as certain non-hematopoietic tissues.² It is activated by the binding of its ligand, stem cell factor (SCF), to the extracellular region, which induces autophosphorylation and the recruitment of signaling molecules to the intracellular region.¹ c-Kit signal transduction activates a variety of signaling pathways, including the PI3K/Akt, RAS/MAPK, and JAK/STAT pathways.^{1,2} The GNNK-isoform of c-Kit undergoes more rapid autophosphorylation and internalization and shows increased MAPK signaling compared with GNNK+ c-Kit, promotes anchorage-independent growth *in vitro*, and induces tumorigenesis in mice.³ Homozygous knockout of *Kit*, the gene encoding c-Kit, is embryonic lethal, with mice developing severe anemia *in utero*.⁴ *KIT* is considered a proto-oncogene and activating mutations have been associated with a variety of cancers.^{1,2} Cayman's c-Kit/CD117 GNNK- Isoform Extracellular Domain (human, recombinant) protein consists of 502 amino acids, has a calculated molecular weight of 56.7 kDa, and a predicted N-terminus of Gln26 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 86 kDa due to glycosylation.

References

1. Roskoski, R., Jr. Src protein-tyrosine kinase structure, mechanism, and small molecule inhibitors. *Pharmacol. Res.* **94**, 9-25 (2015).
2. Edling, C.E. and Hallberg, B. c-Kit—a hematopoietic cell essential receptor tyrosine kinase. *Int. J. Biochem. Cell Biol.* **39(11)**, 1995-1998 (2007).
3. Caruana, G., Cambareri, A.C., and Ashman, L.K. Isoforms of c-KIT differ in activation of signalling pathways and transformation of NIH3T3 fibroblasts. *Oncogene* **18(40)**, 5573-5581 (1999).
4. Russell, E.S. Hereditary anemias of the mouse: A review for geneticists. *Adv. Genet.* **20**, 357-459 (1979).

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