

PRODUCT INFORMATION

K-Ras Isoform B (G12C, C51S, C80L, C118S mutant; human, recombinant)
Item No. 40366

Overview and Properties

Synonyms: c-K-ras(G12C, C51S, C80L, C118S), K-Ras4A(G12C, C51S, C80L, C118S), K-Ras Cys-light, Ki-Ras(G12C, C51S, C80L, C118S), c-Ki-ras(G12C, C51S, C80L, C118S), Kirsten Rat Sarcoma Virus(G12C, C51S, C80L, C118S)

Source: Recombinant human N-terminal His-tagged K-Ras isoform B (G12C, C51S, C80L, C118S) expressed in *E. coli*

Amino Acids: 1-169

Uniprot No.: P01116

Molecular Weight: 22.27 kDa

Storage: -80°C (as supplied)

Stability: ≥1 year

Purity: **batch specific** (≥90% estimated by SDS-PAGE)

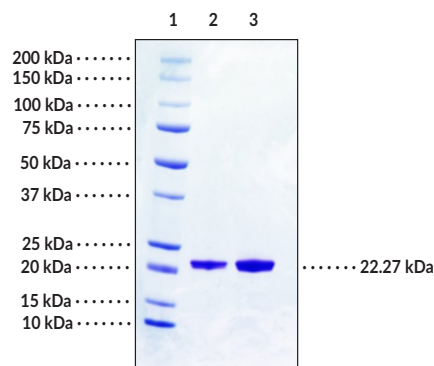
Supplied in: 20 mM HEPES, pH 8.0, with 150 mM sodium chloride, and 1 mM magnesium chloride

Protein

Concentration: **batch specific** mg/ml

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

Image



Lane 1: MW Markers
Lane 2: K-Ras Isoform B (G12C, C51S, C80L, C118S mutant) (2 µg)
Lane 3: K-Ras Isoform B (G12C, C51S, C80L, C118S mutant) (4 µg)

SDS-PAGE Analysis of K-Ras Isoform B (G12C, C51S, C80L, C118S mutant).
This protein has an apparent molecular weight of 22.27 kDa.

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, 12/10/2024

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

K-Ras is a small GTPase and member of the RAS family of GTPases with roles in apoptosis, as well as cell proliferation, survival, and migration.^{1,2} K-Ras is composed of a guanine nucleotide-binding domain containing an active site, an effector-binding domain, and an isoform-specific C-terminal hypervariable region, which varies by four amino acids between isoforms A and B.^{1,3} The active site cycles between GDP-bound inactive and GTP-bound active states and is regulated by its associations with GTPase-activating proteins (GAPs) or guanine nucleotide exchange factors (GEFs).^{3,4} K-Ras is ubiquitously expressed and is tethered to the intracellular side of cell membranes via farnesyl and palmitoyl lipidation.^{1,5} The glycine-to-cysteine substitution at position 12 of mutant K-Ras (K-Ras^{G12C}) is constitutively activating and found in pancreatic, colon, and lung cancers.^{2,6} K-Ras^{G12C} substitutions of serine, leucine, and serine for the native cysteine residues in positions 51, 80, and 118 (C51S, C80L, C118S), respectively, of the guanine nucleotide-binding domain have minimal effects on protein structure, therefore, this tetramutated protein has been used to improve labeling of Cys¹² for protein crystallography and to reduce off-target inhibitor ligation and prevent cysteine-based reactions in kinetic assays.⁷⁻⁹ Cayman's K-Ras Isoform B (G12C, C51S, C80L, C118S mutant; human, recombinant) protein can be used for ELISA, protein crystallography, and Western blot applications.

References

1. Nussinov, R., Tsai, C.-J., Chakrabarti, M., *et al.* A new view of Ras isoforms in cancers. *Cancer Res.* **76**(1), 18-23 (2016).
2. Padavano, J., Henkhaus, R.S., Chen, J., *et al.* Mutant K-RAS promotes invasion and metastasis in pancreatic cancer through GTPase signaling pathways. *Cancer Growth Metastasis* **8**(Suppl 1), 95-113 (2015).
3. Hillig, R.C., Sautier, B., Schroeder, J., *et al.* Discovery of potent SOS1 inhibitors that block RAS activation via disruption of the RAS-SOS1 interaction. *Proc. Natl. Acad. Sci. USA* **116**(7), 2551-2560
4. Mechanism of inhibition by arachidonic acid of the catalytic activity of ras GTPase-activating proteins. *J. Biol. Chem.* **271**(3), 1566-1572 (1996).
5. Salim, A.A., Tan, L., Huang, X.-C., *et al.* Oligomycins as inhibitors of K-Ras plasma membrane localisation. *Org. Biomol. Chem.* **14**(2), 711-715 (2016).
6. Hallin, J., Engstrom, L.D., Hargis, L., *et al.* The KRAS^{G12C} inhibitor MRTX849 provides insight toward therapeutic susceptibility of KRAS-mutant cancers in mouse models and patients. *Cancer Discov.* **10**(1), 54-71 (2020).
7. Ostrem, J.M., Peters, U., Sos, M.L., *et al.* K-Ras(G12C) inhibitors allosterically control GTP affinity and effector interactions. *Nature* **503**(7477), 548-551 (2013).
8. Zhang, Z., Guiley, K.Z., and Shokat, K.M. Chemical acylation of an acquired serine suppresses oncogenic signaling of K-Ras(G12S). *Nat. Chem. Biol.* **18**(11), 1177-1183 (2022).
9. Huynh, M.V., Parsonage, D., Forshaw, T.E., *et al.* Oncogenic KRAS G12C: Kinetic and redox characterization of covalent inhibition. *J. Biol. Chem.* **298**(8), 102186 (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