

PRODUCT INFORMATION



K-Ras(G12D) Isoform A (human, recombinant)

Item No. 40372

Overview and Properties

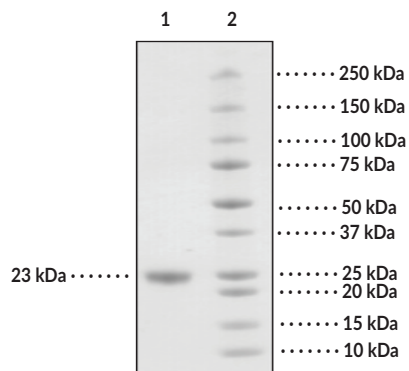
Synonyms: c-K-ras(G12D), c-Ki-ras(G12D), K-Ras4A(G12D), Ki-Ras(G12D), Kirsten Rat Sarcoma Virus(G12D), KRAS(G12D)
Source: Recombinant human N-terminal His-tagged K-Ras(G12D) isoform A expressed in *E. coli*
Amino Acids: 2-186
Uniprot No.: P01116
Molecular Weight: 23 kDa
Storage: -80°C (as supplied)
Stability: ≥6 months
Purity: *batch specific* (≥90% estimated by SDS-PAGE)
Supplied in: 20 mM HEPES, pH 7.4, 150 mM sodium chloride, 10% glycerol, and 1 mM dithiothreitol

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: K-Ras(G12D) Isoform A
Lane 2: MW Markers

SDS-PAGE Analysis of K-Ras(G12D) Isoform A.

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, 03/13/2024

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 (HVR) that differs between K-Ras isoforms A and B due to alternative splicing.^{1,3,4} 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).^{3,5} K-Ras is ubiquitously expressed and is tethered to the intracellular side of cell membranes *via* farnesyl and palmitoyl lipidation and to negatively charged or neutral regions of the membrane *via* a single polybasic region in the HVR.^{1,6,4} K-Ras(G12D), which contains a glycine-to-aspartic acid substitution at position 12, is constitutively active and associated with pancreatic, colon, and lung cancers.⁷ Inhibition of K-Ras(G12D) with the inhibitory peptide KS-58 reduces tumor volume in a PANC-1 pancreatic cancer mouse xenograft model.⁸ Tumor levels of K-Ras(G12D) are increased in patients with lung adenocarcinoma who had never smoked.⁹ Cayman's K-Ras(G12D) Isoform A (human, recombinant) protein consists of 185 amino acids and has a calculated molecular weight of 23 kDa.

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 (2019).
4. Parker, J.A. and Mattos, C. The K-Ras, N-Ras, and H-Ras isoforms: Unique conformational preferences and implications for targeting oncogenic mutants. *Cold Spring Harb. Perspect. Med.* **8(8)**, a031427 (2018).
5. Sermon, B.A., Eccleston, J.F., Skinner, R.H., *et al.* Mechanism of inhibition by arachidonic acid of the catalytic activity of ras GTPase-activating proteins. *J. Biol. Chem.* **271(3)**, 1566-1572 (1996).
6. 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).
7. Swiatnicki, M., Engel, L., Shrestha, R., *et al.* Profiling oncogenic KRAS mutant drugs with a cell-based Lumit p-ERK immunoassay. *SLAS Discov.* **S2472-5552(22)**, (2022).
8. Sakamoto, K., Masutani, T., and Hirokawa, T. Generation of KS-58 as the first K-Ras(G12D)-inhibitory peptide presenting anti-cancer activity in vivo. *Sci. Rep.* **10**, 21671 (2020).
9. Dogan, S., Shen, R., Ang, D.C., *et al.* Molecular epidemiology of EGFR and KRAS mutations in 3,026 lung adenocarcinomas: Higher susceptibility of women to smoking-related KRAS-mutant cancers. *Clin. Cancer Res.* **18(22)**, 6169-6177 (2012).

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