

AbFlex® KRAS antibody (rAb)

Catalog Nos: 91437, 91438

Application(s): ELISA, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Molecular Weight: 24 kDa

Background: AbFlex® antibodies are recombinant antibodies (rAbs) that have been generated using defined DNA sequences to produce highly specific, reproducible antibodies. Each AbFlex antibody contains a 6xHis Tag, a Biotinylation Tag for enzymatic biotin conjugation using the biotin ligase, BirA, and a sortase recognition motif (LPXTG) to attach a variety of labels directly to the antibody including fluorophores, enzymatic substrates (HRP, AP), peptides, drugs as well as solid supports.

AbFlex® KRAS antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

Oncogenic mutations in Ras genes have been found in about 30% of human cancers. This family of genes is involved in cell signaling pathways that control cell growth and death. The human Ras gene family consists of three identified members H-, K- and N-ras encoding proteins of 188-189 amino acids and 21,000 (p21) daltons. Human H- and K-ras are the homologues of v-H- and v-K-ras, sequences. Under normal circumstances the Ras-GTPase activity gradually converts GTP to GDP, but in oncogenic mutations Ras locks the protein into the active GTP conformation. This leads to a constant Ras gene and uncontrollable cell growth. Due to these intracellular relationship Ras antibodies have emerged as a valuable tool for target validation studies for cancer.

Immunogen: This information is proprietary to Active Motif and/or its suppliers.

Buffer: Purified IgG 39mM NaH₂PO₄, 117mM NaCl, 163mM imidazole, 0.035% sodium azide, 30% glycerol

Application Notes:

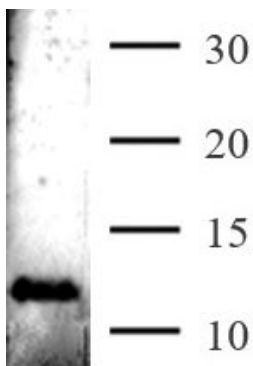
Applications Validated by Active Motif:

WB: 0.5-2 µg/ml dilution

ELISA: Individual optimization may be required.

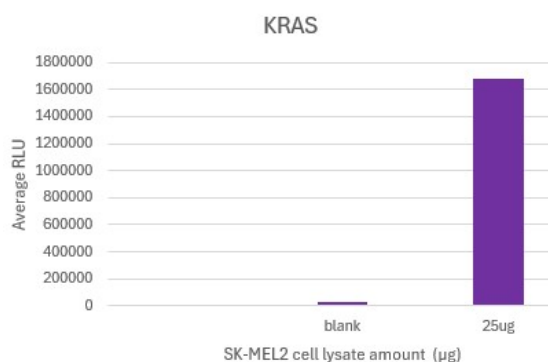
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -80°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



AbFlex® KRAS Antibody (rAb) tested by Western blot.

ASPC-1 WCE whole cell extract (20 µg) was probed with AbFlex® KRAS antibody at 1 mg/ml.



AbFlex® KRAS Antibody tested by bead-based assay multiplex ELISA

Luminex bead-based specificity analysis was used to confirm the specificity of KRAS antibody for recombinant human RAS protein isoforms conjugated to MagPlex Luminex beads. Peptide-bound antibody was detected with anti-mouse IgG-Phycoerythrin and read in a Luminex instrument. (Luminex® is a registered trademark of Luminex Corporation.)