

ARAF Protein

Recombinant human protein expressed in Sf9 cells

Catalog # A24-35G

Lot # S189-1

Product Description

Recombinant human ARAF (209-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_001654](#).

Gene Aliases

PKS2, A-RAF, ARAF1, RAFA1

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

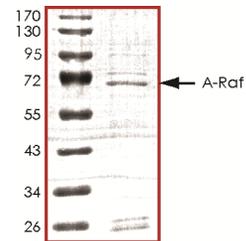
Scientific Background

ARAF belongs to the RAF subfamily of the Ser/Thr protein kinase family, and maybe involved in cell growth and development. ARAF play a critical role in cell growth and development (1).The N-terminal regulatory domain of ARAF interacted with the putative mitochondrial proteins TOM and TIM44. ARAF is expressed predominantly in urogenital tissues (2). The complete coding sequence of the human A-raf-1 oncogene and transforming activity of a human A-raf carrying retrovirus.

References

- Huebner, K. et.al: Actively transcribed genes in the raf oncogene group, located on the X chromosome in mouse and human. Proc. Nat. Acad. Sci. 83: 3934-3938, 1986.
- Lee, J.-E. et.al : The complete sequence and promoter activity of the human A-raf-1 gene (ARAF1). Genomics 20: 43-55, 1994.

Purity



The purity of ARAF was determined to be **>70%** by densitometry. Approx. MW **70 kDa**.

ARAF Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	A24-35G
Specific Lot Number	S189-1
Purity	>70%
Concentration	0.1 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.