

Data Sheet

A-RAF1

Human, recombinant, N-terminal GST tag and C-terminal His tag

Catalog #: 40010

Lot#: 130118-G

Conc.: 0.46 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol and 3 mM DTT.

Stability: >6 months at -80 °C

References:

1. Yokoyama, T. *et al J. Cell Biol.* **177** (5), 781-793. (2007)
2. Ewing, R.M. *et al. Mol. Syst. Biol.* **3**, 89 (2007)

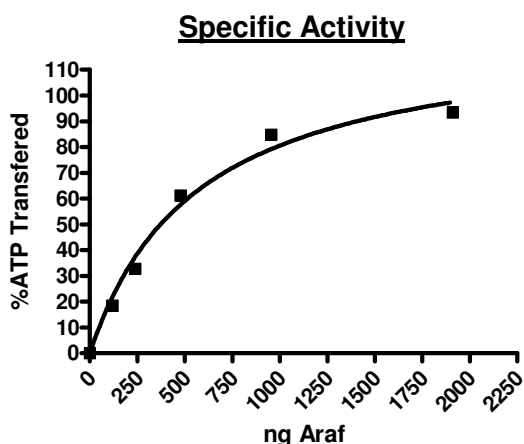
Description: Human v-raf murine sarcoma 3611 viral oncogene homolog, transcript variant 1 (A-RAF1), (GenBank Accession No. NM_001654), amino acids 282-606 (end), containing mutations YY301-302DD, with N-terminal GST tag and C-terminal His tag, MW=63 kDa, expressed in an Baculovirus infected Sf9 cell expression system.

Specific Activity: 40 pmol/min/μg.

Assay conditions: Reaction was done in a kinase buffer (40 mM Tris, 10 mM MgCl₂) containing 5 μg inactive MEK1 (BPS #40075), 1 μM ATP, and 4 – 8 ng/μl aRAF for 30 min at 30°C. The amount of ATP transferred was measured using a Kinase-Glo® Plus Assay Kit (Promega Corporation, Madison, WI).

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

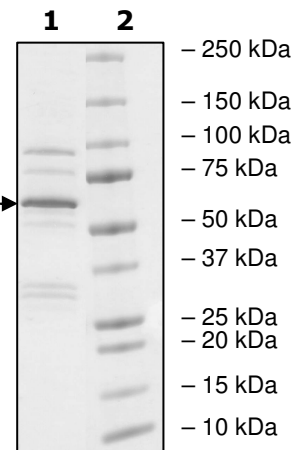


10% SDS-PAGE
Coomassie staining

Lane 1:
2 μg A-RAF1

Lane 2:
Protein Marker

MW: 63 kDa
Purity: ≥62%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.259.5734**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMIIRYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSKDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDPKSLADDKKKVKNLG
YRDSG**DD**WEVPPSEVQLLKRIGTGSFGTVFRGRWHGDVAVKVLKVSQPTAEQAQAF
KNEMQVLRKTRHVNILLFMGMTRPGFAITQWCEGSSLYHHLHVADTRFDMVQLIDV
ARQTAQGMDYLHAKNIIHRDLKSNIFLHEGLTVKIGDFGLATVKTRWSGAQPLEQPS
GSVLWMAAEVIRMQDPNPYSFQSDVYAYGVVLYELMTGSLPYSHIGCRDQIIFMVGR
GYLSPDLSKISSNCPKAMRRLSDCLKFQREERPLFPQILATIELLQRSLPKIERSASE
PSLHRTQADELPACLLSAARLVP**HHHHHH**

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.259.5734**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com