

Data Sheet

ALK

Human, recombinant, N-terminal GST tag
Catalog #: 40214
Lot #: 120919-GC2 **Conc.:** 0.21 mg/ml

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

Stability: >6 months at -80 °C

References:

1. Ren, H., *et al.*, *Cancer Res.* 2012;**72(13)**:3312-3323.
2. Zhu, S., *et al.*, *Cancer Cell* 2012;**21(3)**:362-373.
3. Schonherr, C., *et al.*, *Crit Rev Oncog.* 2012;**17(2)**:123-143.

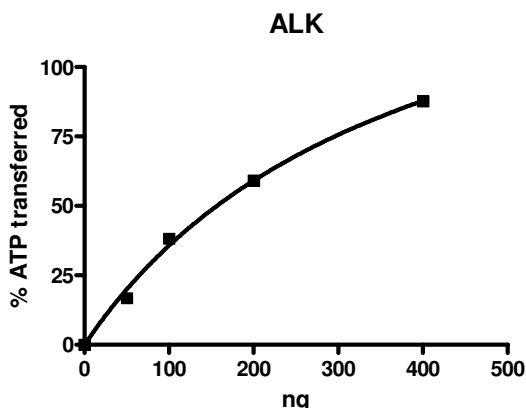
Description: Human anaplastic lymphoma receptor tyrosine kinase (ALK), also known as CD246 and NBLST3, (GenBank Accession No. NM_004304), a.a. 1058-1620 (end) with N-terminal GST tag, MW = 89 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 77 pmole/min/μg.
 Assay conditions: Assay was done in a Kinase buffer using 0.2 mg/ml IGF1Rtide as a substrate with 20 μM ATP at 30°C for 40 min. The amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

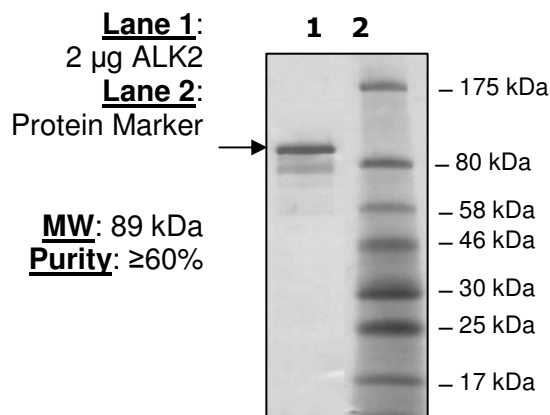
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

Specific Activity



4-20% SDS-PAGE Coomassie staining



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 **1.858.481.8694**
 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
DGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDPLVPRGSPGEFVYRR
KHQELQAMQMELQSPEYKLSKLRSTIMTDYNPNYCFAGKTSSISDLKEVPRKNITLIR
GLGHGAFGEVYEGQVSGMPNDPSPLQVAVKTLPEVCSEQDELDFLMEALIISKFNHQ
NIVRCIGVSLQSLPRFILLELMAGGDLKSFLRETRPRPSQPSSLAMLDLLHVARDIACG
CQYLEENHFHRDIAARNCLLTCPGPGRVAKIGDFGMARDIYRASYYRKGCCAMLPV
KWMPPEAFMEGIFTSKTDTWSEFVLLWEIFSLGYMPYPSKSNQEVLEFVTSGGRMDP
PKNCPGPVYRIMTQCWQHQPEDRPNFAILERIEYCTQDPDVINTALPIEYGPLVEEEE
KVPVRPKDPEGVPPLLVSQQAQKREEERSPAAPPPLPTTSSGKAAKKPTAAEISVRVP
RGPAVEGGHVNMFAFSQSNPPSELHKVHGSRNKPTSLWNPTYGSWFTEKPTKKNPI
AKKEPHDRGNLGLGSCTVPPNVATGRLPGASLLEPSSLTANMKEVPLFRLRHFP
GNVNYGYQQQLPLEAATAPGAGHYEDTILKSKNSMNQPGP

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 **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com