

Data Sheet

ALK1, GST-Tag Active

Human, recombinant, N-terminal GST tag

Catalog #: 40018

Lot#: 180430-A **Conc.:** 0.53 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 0.2mM EDTA, 3mM DTT and 20% glycerol.

Stability: At least 6 months at -80°C . *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

1. Ota, T. et al . J. Cell Physiol. 2002;193(3):299-318.
2. Tsuzuki, T. et al. Am. J. Surg. Pathol. 2004; 28(12):1609-14.

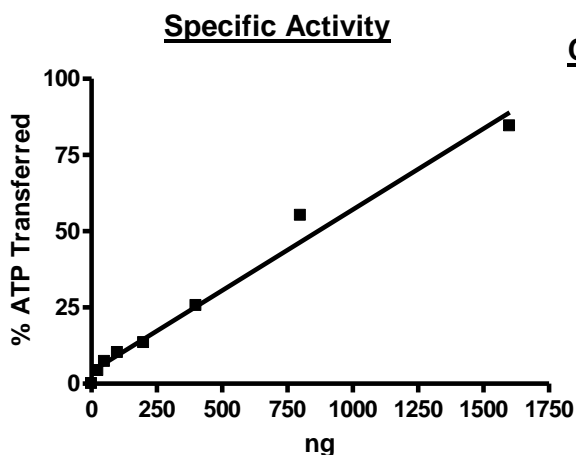
Description: Recombinant human activin receptor-like kinase 1, or ALK1, genbank accession number NM_000020, a.a. 144-503 (end), with N-terminal GST-tag, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 13 pmol/min/ μg

Assay Conditions: Assay was done in Kinase buffer containing 1 mM DTT using Casein substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 45 min. The amount of ATP transferred from autophosphorylation as well as substrate phosphorylation was calculated using Kinase-Glo reagent (Promega).

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

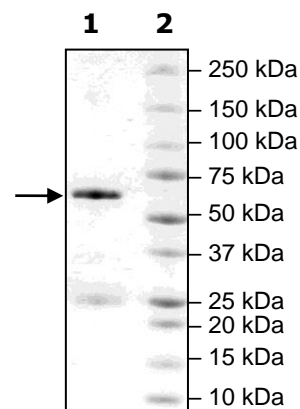
Quality Assurance



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg ALK1
Lane 2:
Protein Marker

MW: 67 kDa
Purity: $\geq 78\%$



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