

Data Sheet

ALK2 (ACVR1), FLAG-Tag, Avi-Tag, Biotin-Labeled Recombinant

Human, Recombinant, N-terminal FLAG-Tag, C-terminal Avi-Tag, Biotin-labeled

Catalog #: 100176

Lot#: 180830-1 **Conc.:** 0.26 mg/ml

Formulated in: 40mM Tris-HCl pH8.0, 110mM NaCl, 2.2mM KCl, 8ng/μL Flag peptide, 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:

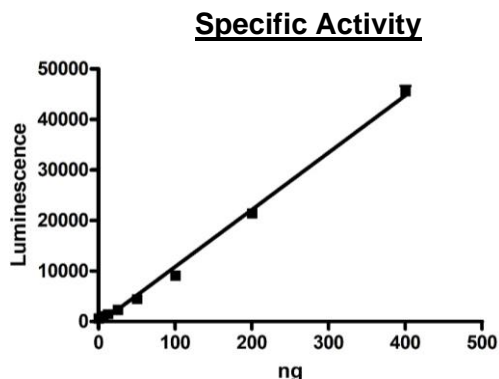
1. Kaplan, F., S., *et al.*, *Hum. Mutat.* 2009;**30**:379-390.
2. Yu, P.B., *et al.*, *Nature Med.* 2008;**14**:1363-1369.

Description: Recombinant human ALK2 a.a. 147-509, (Genbank Accession No. NM_001105.5) with an N-terminal FLAG-tag and C-terminal Avi-Tag™, expressed in Sf9 cells using a baculovirus expression system. MW= 45 kDa.

Assay conditions: Enzyme reactions include 40 mM Tris pH 7.4, 20 mM MgCl₂, 0.1 mg/mL BSA, 10 μM ATP, 0.2 mg/mL casein, 3 mM DTT, and various amounts of ALK2—incubated 45 minutes at 30 °C. Followed by addition of 25 μL of BPS ADP-Glo reagent—incubated 45 minutes at room temperature. Followed by addition of 50 μl of BPS Kinase Detection Reagent—incubated 45 minutes at room temperature. Luminescence was measured.

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

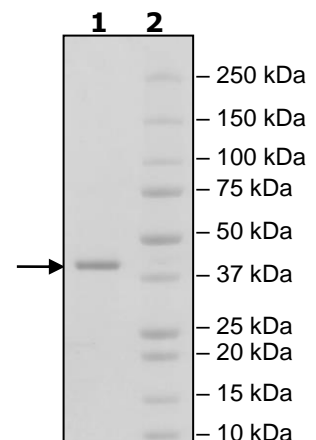
Quality Assurance



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

Lane 1:
3 μg ALK2
Lane 2:
Protein Marker

MW: 45 kDa
Purity: 90%



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com