

Data Sheet

ALK2, FLAG-Tag, Avi-Tag, Biotin-Labeled

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:

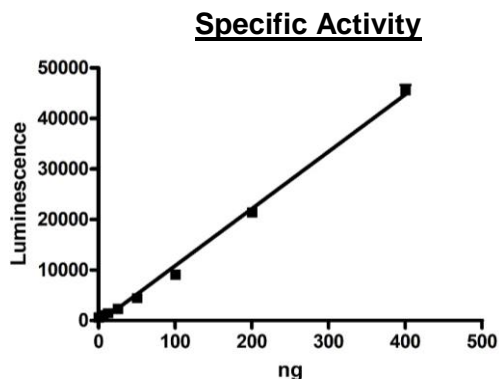
1. Kaplan, F., S., *et al.*, *Hum. Mutat.*
 2009;**30**:379-390.
2. Yu, P.B., *et al.*, *Nature Med.*
 2008;**14**:1363-1369.

Description: Recombinant human ALK2
 a.a. 147-509, (Genbank Accession No.
 NM_1001105) 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.

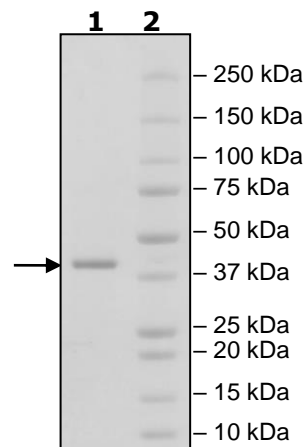
Quality Assurance



**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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com