

Data Sheet

ALK2 GST-Tag

Human, Recombinant, N-terminal GST-tag

Catalog #: 40019

Lot#: 180816

Conc.: 0.63 mg/ml

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

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-end (Genbank Accession No. NM_1001105) with an N-terminal GST-tag, expressed in Sf9 cells using a baculovirus expression system. MW = 67 kDa.

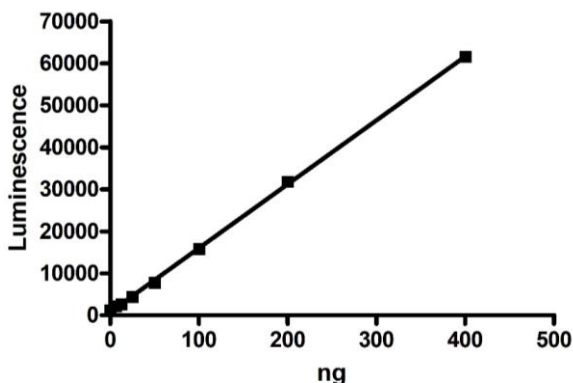
Specific Activity: 40 pmol/min/ μg

Assay Conditions: Enzyme reactions include 40 mM Tris pH 7.4, 20 mM MgCl_2 , 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

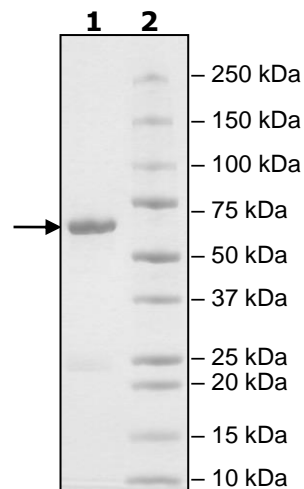
Specific Activity



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg ALK2
Lane 2:
Protein Marker

MW: 67 kDa
Purity: $\geq 90\%$



LY. 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