

Data Sheet

ALK2 (ACVR1), GST-tag Recombinant

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.*

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.

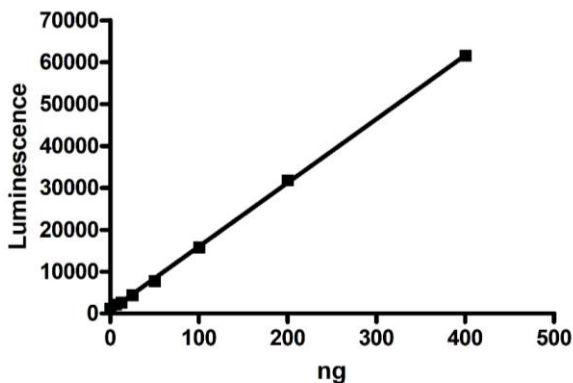
Description: Recombinant human ALK2 a.a. 147-end, Genbank Accession No. NM_001105.5 with an N-terminal GST-tag, expressed in Sf9 cells using a baculovirus expression system. MW = 67 kDa.

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

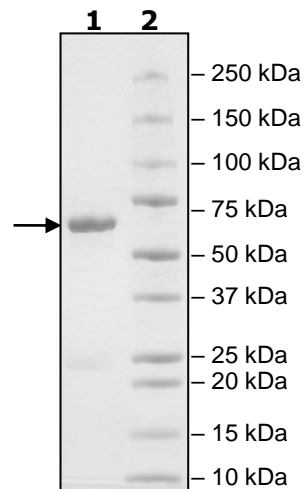


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg ALK2
Lane 2:
Protein Marker

MW: 67 kDa

Purity: $\geq 90\%$



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