

Data Sheet

ULK1, Flag-Tag

Human, recombinant, N-terminal Flag-tag

Catalog #: 40099

Lot #: 150806-B **Conc.:** 0.25 mg/ml

Formulated in: 25 mM Tris-HCl, pH 7.5, 100 mM NaCl, 0.05% Tween-20, 100 µg/ml FLAG peptide, 20% glycerol and 3 mM DTT.

Stability: At least 6 months at -80°C. *Avoid freeze/thaw cycles.*

References:

1. Ganley I.G. et al., J. Biol. Chem. **284 (18)**, 12297-12305 (2009).
2. Chan E.Y. et al., Mol. Cell. Biol. **29 (1)**, 157-171 (2009).

Description:

Human ULK1, also known as ATG1, UNC51, and hATG1, GenBank Accession No. NM_003565, a.a. 2-1050 (end) with N-terminal Flag-tag, expressed in a HEK293 expression system. MW = 114 kDa

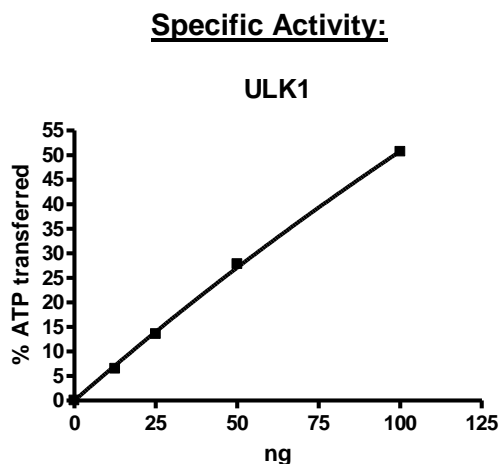
Specific Activity: ≥85 pmol/min/µg

Assay Conditions: Assay was done in a Kinase buffer using MBP (0.2 mg/ml) as a substrate with 20 µM ATP at 30°C for 60 min. The amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

Application:

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

Quality Assurance

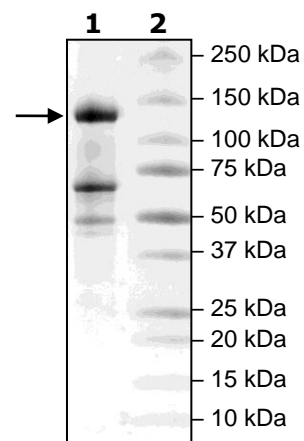


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 µg ULK1

Lane 2:
Protein Marker

MW: 114 kDa
Purity: ≥40%



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