

Data Sheet

VPS34, FLAG-Tag

Human, recombinant, N-terminal FLAG

Catalog #: 40690

Lot#: 180611-2 **Conc.:** 0.13 mg/ml

Formulated in: 40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 µg/mL FLAG peptide, 20% glycerol, 3 mM DTT.

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

References:

1. Yan, Y., *et al.* (2009). *Biochem J.* **417(3):**747-55.
2. Backer, J.M. (2008). *Biochem J.* **410(1):**1-17.

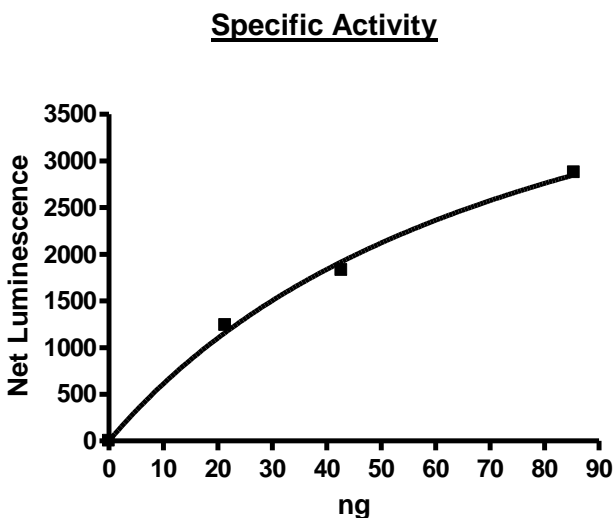
Description: Human VPS34 also known as vacuolar protein sorting 34, PI3-Kinase Type 3, and Phosphatidylinositol 3-Kinase Catalytic Subunit Type 3, GenBank Accession No. NM_002647, a.a. 2-887(end) with N-terminal FLAG-tag, expressed in Sf9 insect cells via a baculovirus expression system. MW=103 kDa.

Specific Activity: 120 pmole/min/µg.

Assay was done in a Kinase buffer containing 2 mM MgCl₂ at 30°C for 20 min. Use BPS 4PI:PS substrate (100 µM) with 100 µM ATP. The amount of ATP transferred was calculated using ADP-Glo reagent.

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

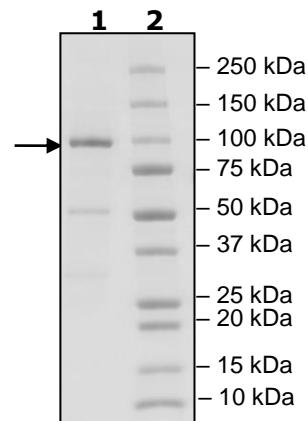
Quality Assurance



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

Lane 1:
3 µg VPS34
Lane 2:
Protein Marker

MW: 103 kDa
Purity: ≥80%



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