

Data Sheet

KAT5 (TIP60), GST-Tag

Human, Recombinant, with N-terminal GST-tag

Catalog #: 79422

Lot #: 180123 **Conc.:** 0.1 mg/ml

Formulated in: 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

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

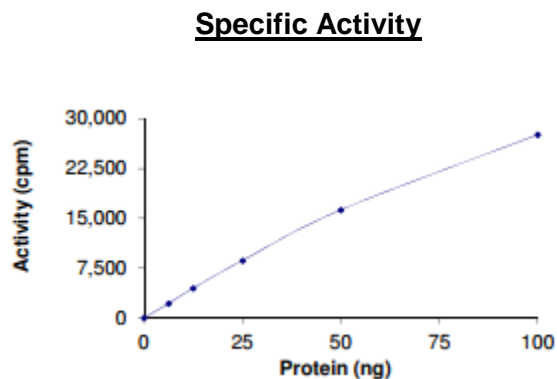
References:

1. Kamine, J. *et.al. Virology.* 1996; **216**: 357-366.
2. Gorrini, C. *et al. Nature.* 2007 **448**:1063-1067.

Description: Full-length recombinant Human KAT5 (TIP60) protein also known as cPLA2 ESA1 HTATIP, HTATIP1, PLIP, and TIP. Genbank Accession NM_006388, with an N-terminal GST tag expressed by baculovirus in Sf9 insect cells. MW = 88kDa.

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

Quality Assurance

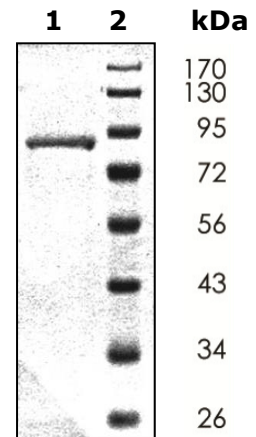


**10% SDS-PAGE
Coomassie staining**

Lane 1:
KAT5 Protein

Lane 2:
Protein Marker

MW: ~88 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