

## Data Sheet

### TYRO3, GST-Tag

Human, Recombinant, N-terminal GST-Tag

**Catalog #:** 40293

**Lot #:** 170427-3

**Conc.:** 0.48 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 1.6mM Glutathione

**Stability:** Store product at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

1. Angelillo-Scherrer, A. *et al. Clin Invest.* 2005; **115**: 237-246.
2. Liao, X., *et al. Mammalian Genome.* 1996; **7**: 395-396.

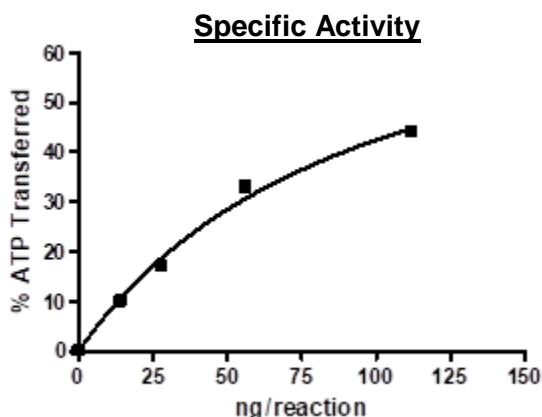
**Description:** Human TYRO3, also known as BYK, DTK, and TYRO3 protein tyrosine kinase, GenBank Accession No. NM\_006293, a.a. 455-890 (end), MW= 74 kDa, expressed in SF9 insect cells.

**Specific Activity:** 140 nmol/min/mg  
 Assay Conditions: Assay was done in Kinase buffer containing 2 mM DTT using Poly-(Glu4:Tyr) substrate (0.2 mg/ml) and 20  $\mu\text{M}$  ATP. Reaction was done at  $30^{\circ}\text{C}$  for 45 min. 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



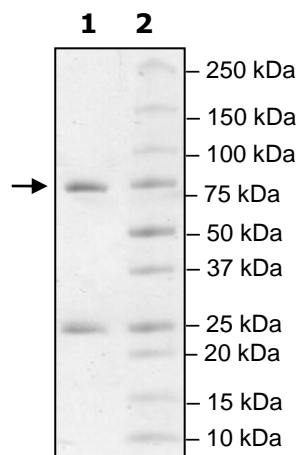
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
TYRO3(455-e)

**Lane 2:**  
Protein Marker

**Purity:**  $\geq 55\%$

**MW:** 74 kDa



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)