

## Data Sheet

### FGFR3, His-Avi-Tag, Biotin-labeled

Human, Recombinant, N-terminal His-Avi-tag,  
Biotin-labeled

**Catalog #:** 100613

**Lot #:** 200130

**Conc:** 0.28 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

#### References:

1. Cappellen, D. *et al.*, *Nature Genetics*. (1999); **23**:18-20.
2. Bellus, G.A. *et al.*, *AM. J. Hum. Genet.* (1995); **56(2)**:368-373.

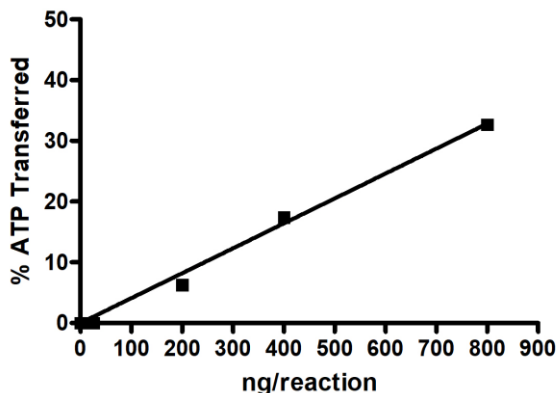
**Description:** Human Fibroblast growth factor receptor 3 (FGFR3), also known as JTK4, CD333, CEK2 and ACH, GenBank Accession No. NM\_000142, a.a. 447-761, with N-terminal His-Avi-tag, expressed in an Sf9 infected baculovirus expression system and enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be  $\geq 90\%$ . MW = 39 kDa. This protein was isolated from a co-expression with PTP1B. Percentage of phosphorylation  $\leq 5\%$ .

**Specific Activity:** 4.3 pmole/min/ $\mu\text{g}$   
**Assay Conditions:** Assay was done in Kinase buffer containing 2 mM DTT using 0.2 mg/mL Poly-(Glu:Tyr) substrate and 10  $\mu\text{M}$  ATP. Reaction was done at  $30^{\circ}\text{C}$  for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

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

## Quality Assurance

### Specific Activity

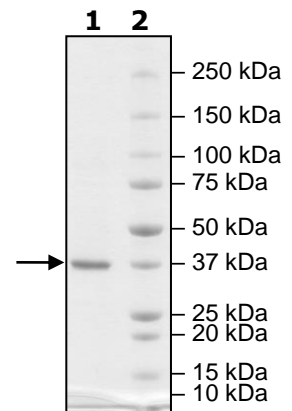


### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2  $\mu\text{g}$  FGFR3  
**Lane 2:**  
Protein Marker

**MW:** 39 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 **1.858.202.1401** 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)