

## Data Sheet

### **FGFR3 (K650E), His-Tag (Dephosphorylated)**

Human, Recombinant, N-terminal His-Tag  
**Catalog #:** 100782  
**Lot#:** 200722                      **Conc.:** 1.96 mg/ml

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

**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, and Hydroxyaryl-Protein Kinase with K650E mutation, GenBank Accession No. NM\_000142, a.a. 447-761, with N-terminal His-tag, co-expressed with PTP1B for dephosphorylation in an Sf9 cell expression system, phosphorylation level <5%.  
MW = 37 kDa.

## **Quality Assurance**

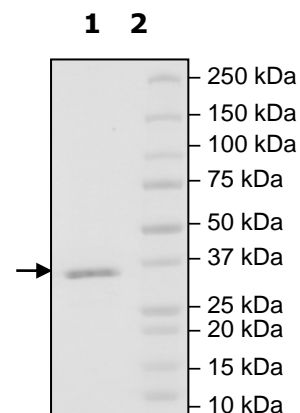
### 4-20% SDS-PAGE Coomassie staining

Lane 1:  
FGFR3-K650E

Lane 2:  
Protein Marker

MW: 37 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 **1.858.202.1401** 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)