

## Data Sheet

### **FGFR3 (V555F), His-Tag (Dephosphorylated)**

Human, Recombinant, N-terminal His-Tag  
**Catalog #:** 100657  
**Lot#:** 200706-G      **Conc.:** 0.51 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:

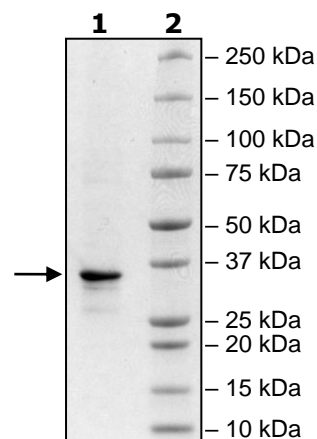
1. Cappellen, D. *et al.*, *Nature Genetics*. (1999); **23**:18-20.
2. Bellus, G.A. *et al.*, *AM. J. Hum. Genet.* (1995); **56(2)**:368-373.

**Description:** Human Fibroblast Growth Factor Receptor 3 (FGFR3), also known as JTK4, and Hydroxyaryl-Protein Kinase with V555F 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

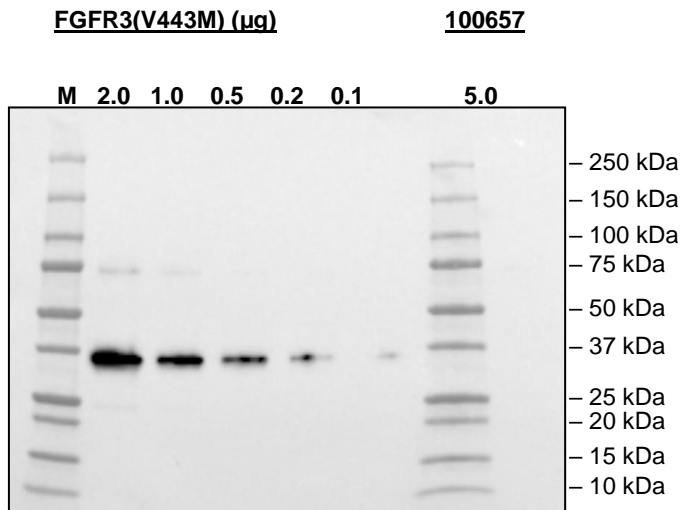
**Lane 1:**  
2 $\mu\text{g}$  FGFR3  
**Lane 2:**  
Protein Marker  
**MW:** 37 kDa  
**Purity:** >90%



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)

**Anti-pTyr Western Blot**



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)