

## Data Sheet

### FGFR3 (V443L), His-Tag

Human, Recombinant, with N-terminal His-tag

**Catalog #:** 100138

**Lot #:** 180808-G

**Conc:** 1.21 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.02% Tween, 20% glycerol, 3mM DTT

**Stability:** At least 6 months 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. Chen L. et al. Front Biosci. 2005; **296(1)10:** 1961-1976.
2. Mhaweche. et al. Eur.J.Surg.Oncol; 2006 ;**32(2):**231-237.

#### **Description:**

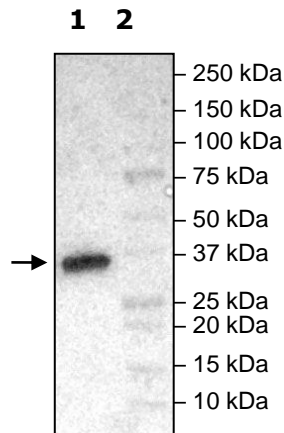
Human FGFR3, GenBank Accession # NM\_000142, with V443L mutation, with N-terminal His-tag, a.a 447-761, MW= 37 kDa, expressed in Sf9 cells via a baculovirus expression system. Phosphorylation is <10%.

### Quality Assurance

#### Western Blot

**Lane 1:**  
FGFR3-V443L

**Lane 2:**  
Protein Marker



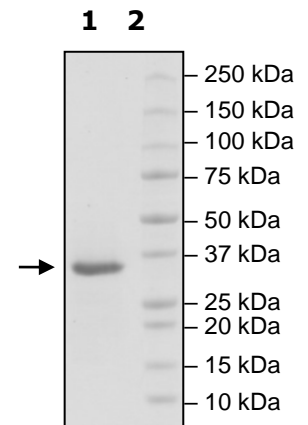
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
FGFR3-V443L

**Lane 2:**  
Protein Marker

**MW:** 37 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 **1.858.829.3082** 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)