

Data Sheet

FGFR3 (L496V), His-Tag

Human, Recombinant, with N-terminal His-tag

Catalog #: 100142

Lot #: 180808-G

Conc: 1.42 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°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. Mhawech. et al. Eur.J.Surg.Oncol; 2006 ;**32(2)**:231-237.

Description:

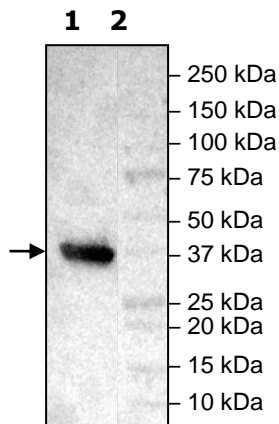
Human FGFR3, GenBank Accession # NM_000142, with L496V mutation, with N-terminal His-tag, a.a 447-761, MW= 37 kDa, expressed in Sf9 cells via a baculovirus expression system. Phosphorylation is <10%.

Quality Assurance

Western Blot

Lane 1:
FGFR3-L496V

Lane 2:
Protein Marker



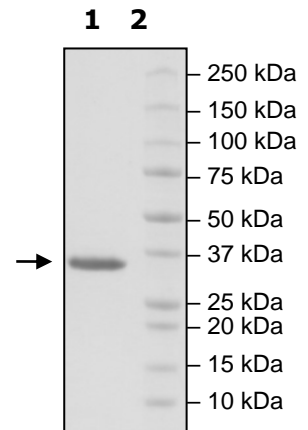
4-20% SDS-PAGE Coomassie staining

Lane 1:
FGFR3-L496V

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

Please visit our website at: www.bpsbioscience.com