

Data Sheet

FGFR3 (447-761) (K650M), His-Tag (Dephosphorylated)

Human, Recombinant, N-terminal His-Tag
Catalog #: 100783
Lot#: 200722 **Conc.:** 1.18 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°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 K650M 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

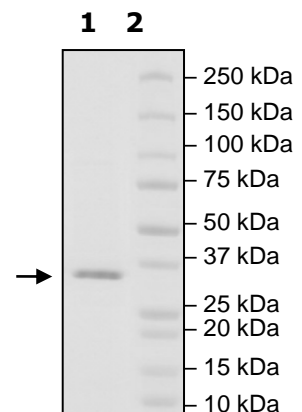
4-20% SDS-PAGE Coomassie staining

Lane 1:
FGFR3-K650M

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: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com