

Data Sheet

FGFR1, Avi-His-Tag, Biotin-labeled (Dephosphorylated) Recombinant

Human, Recombinant, C-terminal Avi-His-tag,
 Biotin-labeled

Catalog #: 100692

Lot #: 200518

Conc: 1.73 mg/ml

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

Stability: At least 6 months at -80°C . Avoid
 freeze/thaw cycles.

References:

1. Weiss, J. *et al.*, *Sci. Trans. Med.* (2010);
2(62):62ra93.
2. Dode, C. *et al.*, *Nature Genetics.*
 (2003);**33:463-465.**

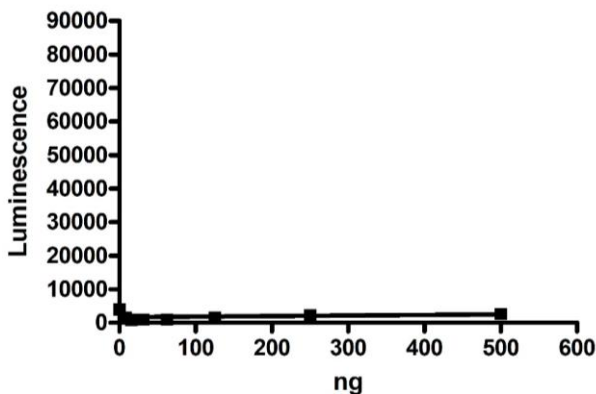
Description: Human Fibroblast growth factor receptor 1 (FGFR1), also known as BFGFR, CD331, CEK and FLT2, GenBank Accession No. NM_023110.3, a.a. 399-822(end), with C-terminal Avi-His-tag, co-expressed with PTP1B for dephosphorylation in an Sf9 cell expression system and enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be $\geq 90\%$ and phosphorylation is $\leq 5\%$. MW=51 kDa.

Assay Conditions: Assay was done according to TYR Phosphorylation Assay Kit (BPS Bioscience #79502) with FGFR enzyme titration range 0 ng/rxn-500 ng/rxn.

Applications: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

FGFR Phosphorylation Detection

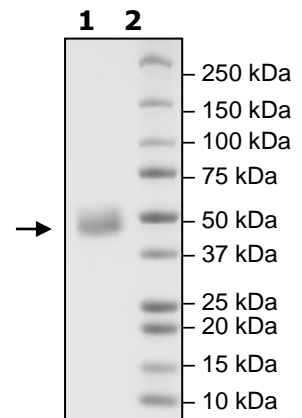


**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
FGFR1
Lane 2:
 Protein Marker

MW: 51 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 **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com