

Data Sheet

FGFR2, DEL (429-430), His-Tag (Dephosphorylated)

Human, Recombinant, C-terminal His-tag

Catalog #: 100356

Lot #: 190502-1

Conc: 0.45 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 100 mM Imidazole, 3 mM DTT and 20% glycerol.

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

References:

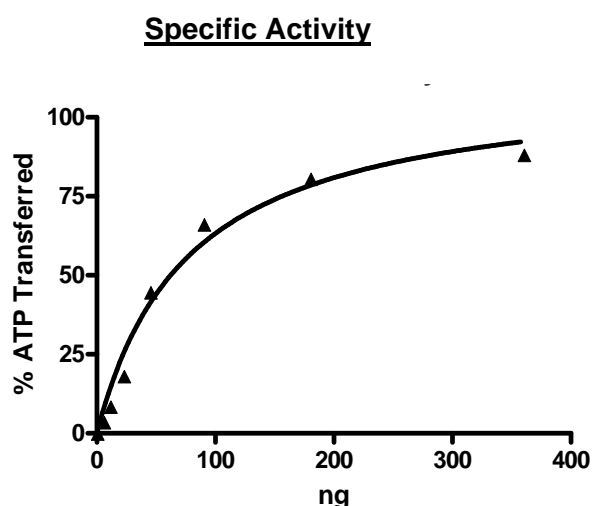
- Holtrich, U. *et al.*, *Proc. Nat. Acad. Sci.* (1991); **88**:10411-10415.
- Roidl, A. *et al.*, *Clin Cancer Res.* (2009 Mar 15); **15**(6):2058-66.

Description: Human FGFR2, also known as Fibroblast growth factor receptor 2, Keratinocyte growth factor receptor, BBDS, BEK, BFR-1 and CD332, GenBank Accession No. NM_000141, a.a. 400-821(end), with deletion at 429-430, with C-terminal His-tag, expressed in an Sf9 infected baculovirus expression system. MW = 52 kDa. This protein was isolated from a co-expression with PTP1B. Percentage of phosphorylation $\leq 10\%$.

Specific Activity: 160 pmole/min/ μg
Assay Conditions: Assay was done according to BPS Catalog#79502 with a FGFR2 titration range of 0 ng/rxn-500 ng/rxn.

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

Quality Assurance



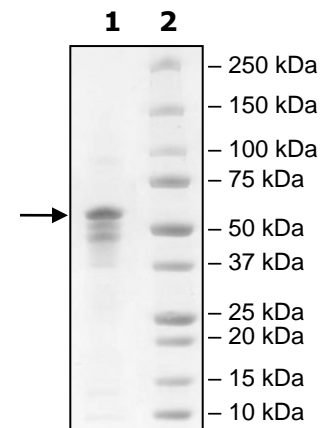
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg FGFR2

Lane 2:
Protein Marker

MW: 52 kDa

Purity: $\geq 70\%$



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

Please visit our website at: www.bpsbioscience.com