

Data Sheet

FGFR2, GST-Tag

Human, Recombinant, N-terminal GST-tag
Catalog #: 40211
Lot #: 171109-GA **Conc:** 0.46 mg/ml

Formulated in: 25mM Tris-HCl, pH 8.0, 138 mM NaCl, 0.05% Tween-20, 50% glycerol and 3mM DTT.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

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.

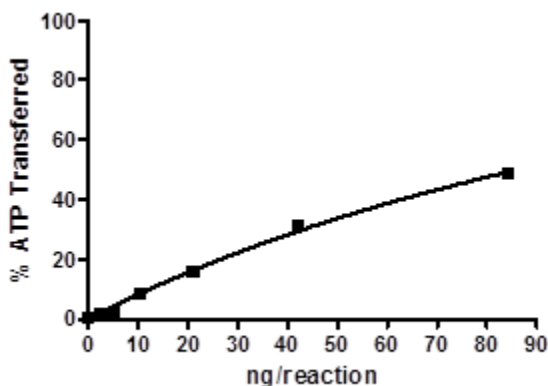
Description: Human FGFR2 (GenBankAccession No. NM_000141), (a.a. 399-end) with N-terminal GST tag, MW= 74 kDa, expressed in baculovirus expression system.

Specific Activity: $\geq 150 \text{ pmol/min}/\mu\text{g}$.
Assay Condition: Assay was done in Kinase buffer containing 2 mM DTT using Poly-(Glu:Tyr) substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

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

Quality Assurance

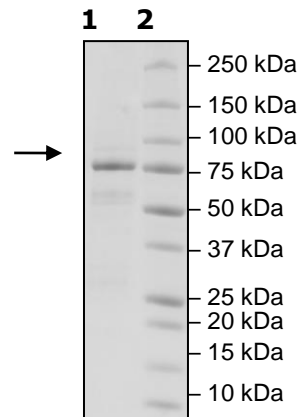
Specific Activity



4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg FGFR2
Lane 2:
Protein Marker

MW: 74 kDa
Purity: $\geq 73\%$



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