

Data Sheet

FGR

Human, recombinant, N-terminal GST tag

Catalog #: 40422

Lot #: 110419

Conc.: 0.1 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.1 mM EDTA, 0.1 mM EGTA, 0.25 mM DTT, 0.1 mM PMSF, 25% glycerol.

Stability: >12 months at -80°C

References:

1. Cheah, M.S., *et al. Nature*. **319(6050)**: 238-240 (1986).
2. Willman, C.L., *et al. Proc. Natl. Acad. Sci. USA* **84(13)**: 4480-4484 (1987).

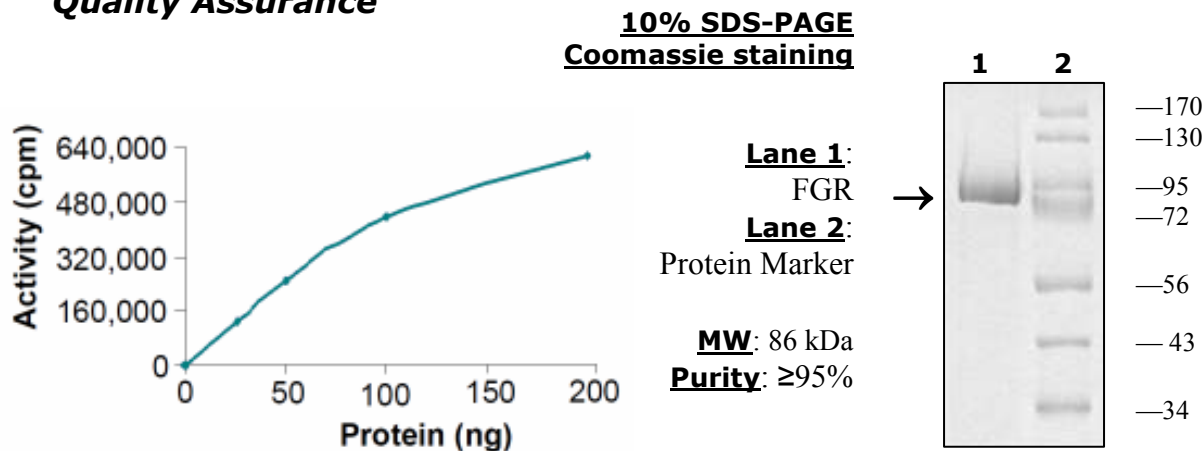
Description: Human FGR, also known as SRC2, GenBank Accession No. NM_005248, full length, with N-terminal GST tag, MW ~86 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 289 pmol/min/μg

Assay Buffer: 5 mM MOPS, 2.5 mM β-glycerophosphate, 5 mM MgCl₂, 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 2 mM ATP. Incubate FGR with 0.5 mg/ml Poly (Glu:Tyr, 4:1) peptide substrate and [³³P]-ATP at 30°C for 15 minutes, then spot reaction on phosphocellulose paper, fix in 1% phosphoric acid, and assay with a scintillation counter.

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

Quality Assurance



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