

Data Sheet

IRR, GST-Tag

Human, Recombinant, N-terminal GST-Tag

Catalog #: 40242

Lot #: 181210

Conc.: 0.1 mg/ml

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

Stability: At least 6 months at -70°C . Avoid freeze/thaw cycles.

References:

1. Zhang, B. *et al.*, *J. Biol. Chem.* 1992; **267(26)**: 18320-8.
2. Elmlinger, M.W. *et al.*, *Regul. Pept.* 1999. **84(1-3)**: 37-42.

Description:

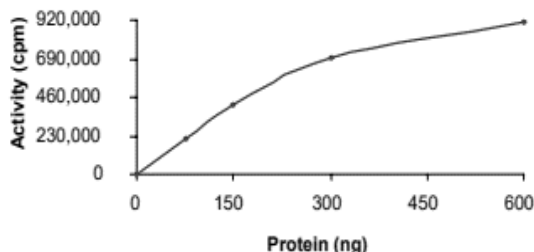
Human IRR protein, also known as Insulin Receptor-related Receptor, and INSR, GenBank Accession No. NM_014215, a.a. 945-end, with N-terminal GST-Tag, expressed in an Sf9 insect cell using a baculovirus expression system. MW = 65 kDa.

Specific Activity: 111 pmol/min/ μg

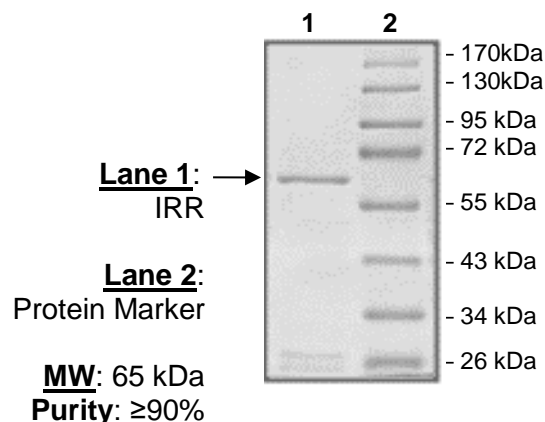
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling

Quality Assurance

Specific Activity



4-20% SDS-PAGE Coomassie Staining



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