

Data Sheet

ROR2, GST-tag

Human, recombinant, N-terminal GST tag
Catalog #: 40296
Lot #: 150723 **Conc.:** 0.10 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 0.1 mM PMSF, and 25% glycerol.

Stability: At least 6 months at -80°C. *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

1. DeChiara, T.M. *et al.*, *Nature Genet.* 2000; **24**: 271-274.
2. Nomi, M. *et al.*, *Molec. Cell Biol.* 2001; **21**: 8239-8335.

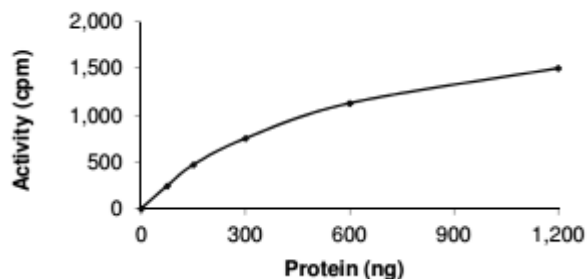
Description: Human ROR2 (GenBank Accession No. NM_004560), a.a. 427-end, with N-terminal GST tag, MW 86 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 0.2 pmol/min/μg

Assay conditions: 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, and 0.25 mg/ml Myelin Basic Protein substrate. Incubate with 0.5 μg ROR1 and 1 μCi [³³P]-ATP at 30°C for 15 minutes, then spot 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.

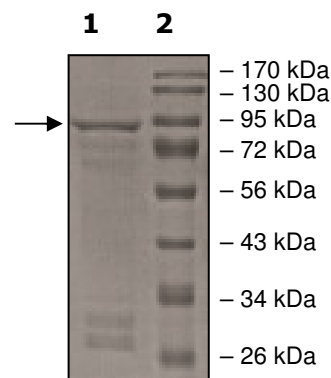
Quality Assurance



**10% SDS-PAGE
 Coomassie staining**

Lane 1: ROR2
Lane 2: Protein Marker

MW: 86 kDa
Purity: ≥80%



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