

Data Sheet

EGFR (D770GY), GST-Tag

Human, Recombinant, with N-terminal
 GST-tag

Catalog #: 79442

Lot #: 180202

Conc.: 0.1 mg/ml

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

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

References:

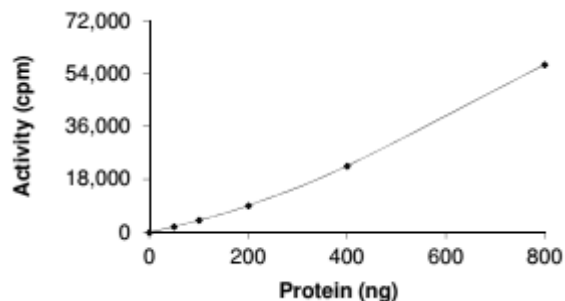
1. Wang, K., *et al.*, *J Biol. Chem*
 2004;**279**:53848-53856.
2. Kobayashi, S., *et al.*, *New Eng. J*
Med. 2005;**325**:786-792.

Description: Recombinant Human
 EGFR also known as ERBB, mENA,
 ERBB1, and HER1 with D770GY
 mutation. Genbank Accession
 NM_005228, a.a. 695 end, with an N-
 terminal GST-tag, expressed in Sf9 cells
 via a baculovirus expression system. MW
 = 89 kDa.

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

Quality Assurance

Specific Activity

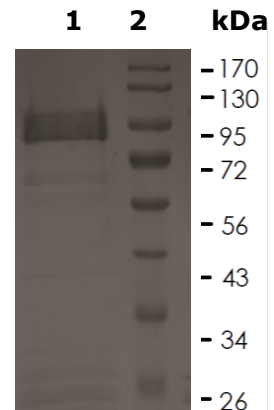


10% SDS-PAGE
Coomassie staining

Lane 1:
 EGFR (D770GY), GST-
 Tag

Lane 2:
 Protein Marker

MW: 89 kDa
Purity: $\geq 80\%$



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