

Data Sheet

EGFR (C775S T790M L858R), GST-Tag

Human, Recombinant, N-terminal
 GST-tag
Catalog #: 79428
Lot #: 180202 **Conc.:** 0.1 mg/ml

Description: Recombinant Human EGFR also known as ERBB, mENA, ERBB1, and HER1 with C775S, T790M, and L858R mutations. Genbank Accession NM_005228, a.a. 668-end, with an N-terminal GST-tag, expressed in Sf9 cells via a baculovirus expression system. MW = 100 kDa.

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.

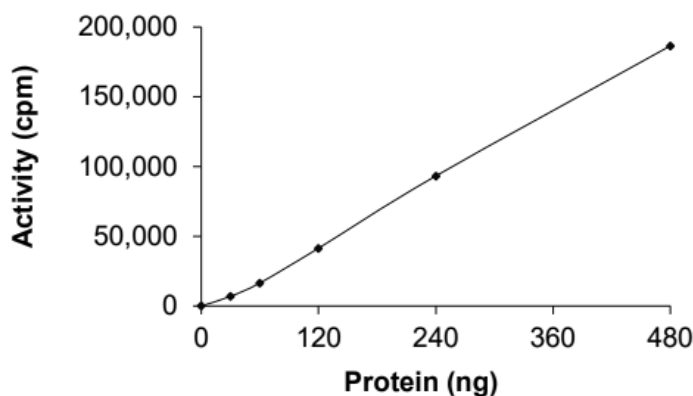
Applications: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

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.

Quality Assurance

Specific Activity



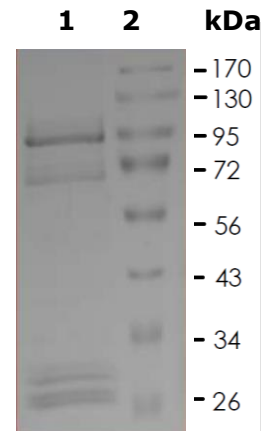
10% SDS-PAGE
Coomassie staining

Lane 1:
 EGFR (C775S T790M L858R) GST-Tag

Lane 2:
 Protein Marker

MW: 100 kDa

Purity: ≥70%



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