

Data Sheet

EGFR (C775S T790M L858R), DEL 746-750, GST-Tag

Human, Recombinant, with N-terminal
GST-tag
Catalog #: 79431
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:

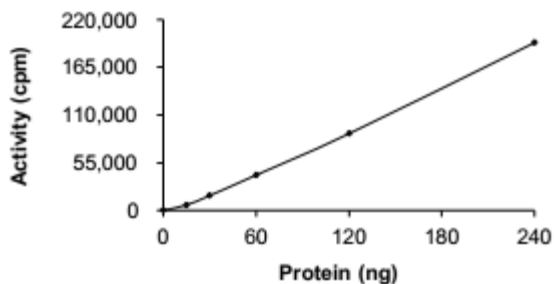
1. Wang, K., *et al.*, *J Biol. Chem*
2004;**279**:53848-53856.
2. Kobayashi, S., *et al.*, *New Eng. J*
Med. 2005;**325**:786-792.

Description: Recombinant Human
EGFR also known as ERBB, mENA,
ERBB1, and HER1 with C797S, T790M,
and L858R mutations and 746-750
deletion. 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.

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

Quality Assurance

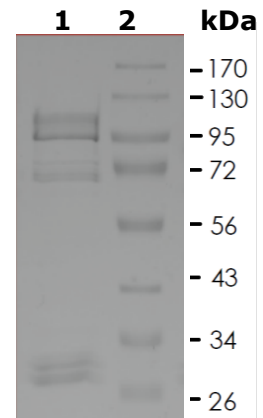
Specific Activity



10% SDS-PAGE Coomassie staining

Lane 1:
EGFR (C775S T790M
L858R), DEL 746-750,
GST-Tag
Lane 2:
Protein Marker

MW: 100 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