

Data Sheet

EGFR (C797S), DEL 746-750, GST-Tag

Human, Recombinant, with N-terminal
GST-tag
Catalog #: 79433
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:

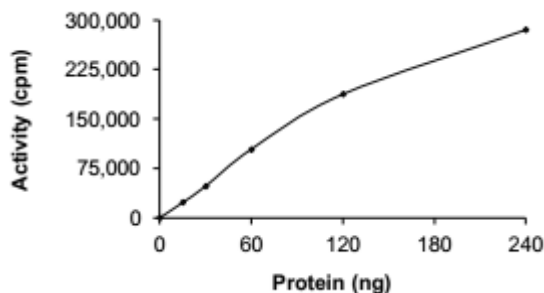
1. Wang, K., *et al.*, *J Biol. Chem*
2004;**279**:53848-53856.
2. Kobayashi, S., *et al.*, *New Eng. J*
Med. 2005;**325**:786-792.

Description: Recombinant Human
EGFR also known as ERBB, mENA,
ERBB1, and HER1 with C797S 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

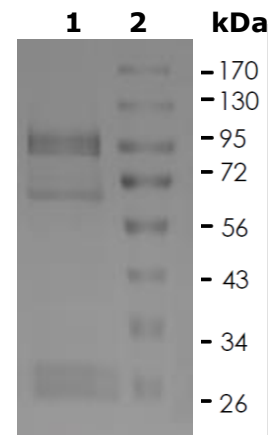


10% SDS-PAGE **Coomassie staining**

Lane 1:
EGFR (C797S), DEL
746-750, GST-Tag

Lane 2:
Protein Marker

MW: 100 kDa
Purity: $\geq 75\%$



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