

PRODUCT DATA SHEET



PEPTIDE NAME: EGFR(988-998), PTP1B substrate

SEQUENCE: Asp-Ala-Asp-Glu-pTyr-Leu-Ile-Pro-Gln-Gln-Gly

SEQUENCE DESCRIPTION: Sequence is from an autophosphorylation site (Tyr-992) of epidermal growth factor receptor (EGFR).

CATALOG NO.: P-323

LOT NO.: temp

MOLECULAR WEIGHT: 1330.3

PURITY: >97% by HPLC analysis.

AMINO ACID ANALYSIS AND IDENTITY: Confirms.

PEPTIDE CONTENT: 93%

SUPPLIED AS: Trifluoroacetate salt; 1 mg net peptide/vial.

PHYSICAL APPEARANCE: Lyophilized solid.

SOLUBILITY: H₂O

STORAGE: -20°C

USAGE: Excellent substrate for mammalian PTP1B^{1,2,3} (BIOMOL cat.# SE-332) with $K_m=3.9$ μ M and *Yersinia* PTPase (BIOMOL cat.# SE-330). May be used in non-radioactive phosphatase assays incorporating the BIOMOL Green™ reagent (BIOMOL cat.# AK-111)^{4,5}.

OTHER: See attached Guide to Handling and Storing Peptides.

LITERATURE REFERENCES:

M. Sarmiento *et al.* *J. Biol. Chem.* 1998 **273** 26368

Z.-Y. Zhang *et al.* *Anal. Biochem.* 1993 **211** 7

Z.-Y. Zhang *et al.* *Proc. Natl. Acad. Sci. USA* 1993 **90** 4446

B. Martin *et al.* *J. Biol. Chem.* 1985 **260** 14932

K.W. Harder *et al.* *Biochem. J.* 1994 **298** 395