

Data Sheet

Protein Tyrosine Kinase Substrate (poly-Glu,Tyr 4:1)

Ready to use PTK substrate (50X)

Catalog #: 40217 **Lot #:** 191001

Size: 100 µl @ 10 mg/ml (50X)

Formulation: Proprietary buffer for peptide reconstitution

Stability: At least 6 months at -20°C.

References:

1. Varkondi E. *et al.*, *J of Receptors and Signal Transduction*, 2005; **25(1)**; 45-56.

Description:

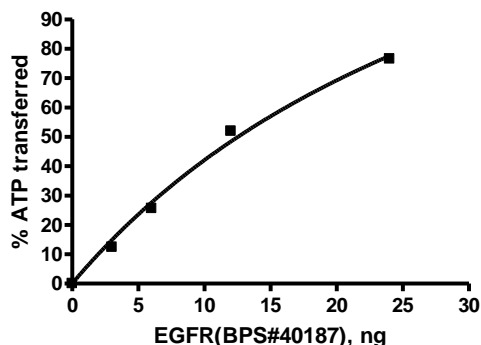
Poly-(Glu,Tyr 4:1)peptide is reconstituted in BPS proprietary buffer. 0.2 mg/ml at final concentration is enough for a majority of PTKs.

Application:

Useful as a substrate for protein tyrosine kinase inhibitor screening.

Quality Assurance

PTK Substrate, poly (Glu, Tyr 4:1)



Enzyme reaction was performed at 30°C for 40 min utilizing 20 µM ATP and PTK substrate (0.2 mg/ml at final). Luminescence signal was generated using Kinase-Glo max reagent (Promega).

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