

## Product Information

<b>Description:</b>	Poly-(Glu,Tyr 4:1) peptide is reconstituted in BPS Bioscience proprietary buffer. 0.20 mg/ml at final concentration is enough for a majority of PTKs.
<b>Concentration:</b>	10 mg/ml
<b>Formulation:</b>	Proprietary buffer for peptide reconstitution
<b>Stability:</b>	At least 6 months at -20°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-20°C
<b>Assay Conditions:</b>	Enzyme reaction was performed at 30°C for 40 min utilizing 20 µM ATP and PTK substrate (0.20 mg/ml at final). Luminescence signal was generated using Kinase-Glo max reagent (Promega).
<b>Applications:</b>	Useful as a substrate for protein tyrosine kinase inhibitor screening.

## Quality Control Data

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

