

## **Product Data Sheet**

Product Name: Jak3tide

Catalog Number: AS-62537 (1 mg) Lot Number: See label on vial

Sequence: H-Gly-Gly-Glu-Glu-Glu-Tyr-Phe-Glu-Leu-Val-Lys-Lys-Lys-OH (3-

letter code)

GGEEEEYFELVKKKK (1-letter code)

Molecular Weight: 1813.0

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Jak3tide is freely soluble in water.

Storage: Jak3tide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at – 20°C or lower. Reconstituted peptide can be aliquoted and stored at –20°C or lower.

Description: This peptide is a substrate for Jak3. It may be used used in kinase assays. Jak3tide contains the phosphorylation site at Tyr7. Misra, U. et al. *J. Biol. Chem.* **281**, 13694 (2006).

For Research Use Only