

Product Data Sheet

| Abltide | |
|---|--|
| AS-61749 (1 mg) | Lot Number: See label on vial |
| Lys-Lys-Gly-Glu-Ala-Ile-Tyr-Ala-Ala-Pro-Phe-Ala-NH2 (3-letter code) KKGEAIYAAPFA-NH2 (1-letter code) | |
| 1264.5 | |
| >95% | |
| Lyophilized white powder | |
| | AS-61749 (1 mg) Lys-Lys-Gly-Glu-Ala-Ile-Tyr-Ala-Ala KKGEAIYAAPFA-NH2 (1-letter cod 1264.5 >95% |

Peptide Reconstitution: Abltide peptide is freely soluble in water.

Storage: Abltide peptide 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: Abltide is a peptide substrate for Abl Kinase (Abl protein tyrosine kinase), a partner in the gag-Abl fusion protein of the Abelson murine leukemia virus. Used in Western blot and kinase assay. Ref: Reddy, P. et al. *Proc. Natl. Acad. Sci. USA.* **80**, 3623 (1983).

Related Products:

| Name | Cat # | Size |
|---|----------|------|
| Abltide, TAMRA-labeled, Abs/Em 541/568 nm | AS-61750 | 1 mg |
| 5-TAMRA-KKGEAIYAAPFA-NH2 | | |

For Research Use Only