

Product Data Sheet

Product Name: Transportan

Catalog Number: AS-61256 (1 mg) Lot Number: See label on vial

Sequence: Gly-Trp-Thr-Leu-Asn-Ser-Ala-Gly-Tyr-Leu-Leu-Gly-Lys-Ile-Asn-Leu-Lys-

Ala-Leu-Ala-Leu-Ala-Lys-Lys-Ile-Leu (3-letter code) GWTLNSAGYLLGKINLKALAALAKKIL (1-letter code)

Molecular Weight: 2841.5

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Transportan peptide is freely soluble in water.

Storage: Transportan 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.

For Research Use Only