



Product Data Sheet

Product Name: TET 830 modified/T-helper epitope from tetanus toxoid

Catalog Number: AS-60622 (1 mg) **Lot Number:** See label on vial

Sequence: H-Ala-Gln-Tyr-Ile-Lys-Ala-Asn-Ser-Lys-Phe-Ile-Gly-Ile-Thr-Glu-Leu-OH
(3-letter code)
AQYIKANSKFIGITEL (1-letter code)

Molecular Weight: 1797.2

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: TET peptide is freely soluble in DMSO.

Storage: TET 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: TET 830 modified/T-helper epitope from tetanus toxoid is a universal human tetanus toxin T cell epitope. It induces T-cell activation and is used as a helper peptide in vaccinations. Ref: Valmori, D. et al. J. Immunol. 149, 717 (1992); Slingluff, C. Jr. et al. Clin. Cancer Res. 7, 3012 (2001); Slingluff, C. Jr. et al. J. Clin. Oncol. 21, 4016 (2003); Demotz, S. et al. J. Immunol. 142, 394 (1989).

For Research Use Only