

## **Product Data Sheet**

Product Name: [Asp370]-Tyrosinase (368-376)

Catalog Number: AS-61456 (1 mg) Lot Number: See label on vial

Sequence: H-Tyr-Met-Asp-Gly-Thr-Met-Ser-Gln-Val-OH (3-letter code)

YMDGTMSQV (1-letter code)

Molecular Weight: 1031.2

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: [Asp370]-Tyrosinase peptide is freely soluble in DMSO.

Storage: [Asp370]-Tyrosinase 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: This tumor antigenic peptide presented by HLA-A2 antigen to CTLs, is a tyrosinase fragment. Tyrosinase is a membrane-bound protein involved in the melanin synthesis pathway that is expressed by virtually all primary melanoma lesions and by most of metastatic lesions. Ref: Chapiro, J. et al. *J. Immunol.* **176**, 1053 (2006); Paczesny, S. et al. *JEM* **199**, 1503 (2004); Lee, P. et al *J. Clin. Oncol.* **19**, 3836 (2001).

For Research Use Only

For Research Use Only