

Product Data Sheet

Product Name: α-Mating Factor Pheromone, yeast

Catalog Number: AS-60221-5 (5 mg) Lot Number: See label on vial

Sequence: H-Trp-His-Trp-Leu-Gln-Leu-Lys-Pro-Gly-Gln-Pro-Met-Tyr-OH (3-letter

sequence)

WHWLQLKPGQPMY (1-letter code)

Molecular Weight: 1684.0 % Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Use distilled or higher quality water. Add water directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5 mg/mL to 1mg/mL or less. Gently vortex to mix. For peptides that have poor solubility in the suggested solvent, brief sonication may increase solubility in some cases.

Storage: Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at – 20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.

Description: This tridecapeptide, an α-factor pheromone of *Saccharomyces cerevisiae*, induces conjugation in yeast by binding to Ste2p. Its cognate GPCR activates a G protein signal pathway that highly conserves with mammalian signaling pathways. Ref: Lee, BK. et al. *J. Biol. Chem.* **276**, 37950 (2001); Mayrhofer, S. and S. Pöggeler *Eukaryot. Cell* **4**, 661 (2005).

For Research Use Only