

Product Data Sheet

Product Name: Orexin B, mouse

Catalog Number: AS-65574 (1 mg) Lot Number: See label on vial

Sequence: Arg-Pro-Gly-Pro-Pro-Gly-Leu-Gln-Gly-Arg-Leu-Gln-Arg-Leu-Leu-Gln-Ala-

Asn-Gly-Asn-His-Ala-Ala-Gly-Ile-Leu-Thr-Met-NH2

RPGPPGLQGRLQRLLQANGNHAAGILTM-NH2

Molecular Weight: 2937.4

Peptide Purity: Peak Area by HPLC ≥95%

Storage: This 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: Orexins are synthesized by posterolateral neurons in hypothalamus. They

bind to orexin-1 (OxR1) and orexin 2 (OxR2) receptors. Orexin A can bind to both OxR1 and OxR2 while Orexin B can associate with OxR2 only. Both Orexins are involved in the regulation of arousal and sleep states, energy

homeostasis, and autonomic central control.

Related Products: Orexin B, human, Cat# AS-65573

Orexin A, Cat# AS-24470

References:

1. Martynska L., et.al., *Neuroendocr Lett* **4(26)**, 289-292 (2005)

2. Ebrahim I.O., et.al., J of Royl Soc of Med 95, 227-230 (2002)

3. Chiou L-C., et.al., Cur Pharm Des 16(28), 3089-3100 (2010)

For Research Use Only