



## Product Data Sheet

**Product Name:** Anti-BetaGamma (MPS-Phosducin-like protein C terminus)

**Catalog Number:** AS-24177 (0.5 mg)      Lot Number: See label on vial  
AS-24178 (1 mg)

**Sequence:** H-Ala-Ala-Val-Ala-Leu-Leu-Pro-Ala-Val-Leu-Leu-Ala-Leu-Leu-Ala-Val-Thr-Asp-Gln-Leu-Gly-Glu-Asp-Phe-Phe-Ala-Val-Asp-Leu-Glu-Ala-Phe-Leu-Gln-Glu-Phe-Gly-Leu-Leu-Pro-Glu-Lys-Glu-OH (3-letter code)  
AAVALLPAVLLALLAVTDQLGEDFFAVDLEAFLQEFGLLPEKE  
(1-letter code)

**Molecular Weight:** 4601.4

**Peptide Purity:** >95%

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:** Initially, try to dissolve a small amount of Anti-BetaGamma peptide in water. If peptide does not dissolve, Anti-BetaGamma peptide is freely soluble in 1% NH<sub>4</sub>OH.

**Storage:** Anti-BetaGamma 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.

**Reference:** Orr, A. et al. *J. Biol. Chem.* **277**, 20453 (2002).

**Additional Information:** Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

The anti-βγ peptide was obtained from AnaSpec, Inc (San Jose, CA). PC12-H<sub>3</sub> and PC12-H<sub>3</sub>/β-ARK1 cells cultured in 12-well plates and differentiated with NGF (100 ng/ml) for 5 to 7 days were incubated with drugs for 5 min. When CBP or the anti-βγ peptide was used, cells were incubated with these compounds for 5 min before incubation with imetit. For pertussis toxin (PTX) treatment, cells were incubated for 24 h with PTX (200 ng/ml) before assay. Dopamine exocytosis was elicited by incubating samples for 5 min with K<sup>+</sup> 100 mM (osmolarity was maintained constant by adjusting the NaCl concentration), phorbol 12-myristate 13-acetate (PMA; 300 nM), or for 20 min with forskolin (10 μM)- [Morrey, C. et al. \*J Pharmacol Exp Ther\* \*\*326\*\*, 871 \(2008\).](#)

### Published Citations:

- Orr, A. et al. *J Biol Chem.* **277**, 20453 (2002).  
Talaia, C. et al. *Neurochem. Intern.* **47**, 418 (2005).  
Morrey, C. et al. *J Pharmacol Exp Ther* **326**, 871 (2008).

*For Research Use Only*