

## **Product Data Sheet**

Product Name: W Peptide, WKYMVm-NH2

Catalog Number: AS-27069 (1 mg) Lot Number: See label on vial

Sequence: Trp-Lys-Tyr-Met-Val-D-Met-NH2 (3-letter code)

WKYMVm-NH2 (1-letter code)

Molecular Weight: 856.1

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: W Peptide is freely soluble in water.

Storage: W 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 peptide, containing the consensus sequence XKYX(P/V)M, is found to stimulate phospholipase C (PLC)-mediated formation of InoPs in certain cell lines and human neutrophils. WKYMVm-NH2 may have the ability to activate the microbicidal functions of human neutrophils. Ref: Baek, S. et al. *J. Biol. Chem.* **271**, 8170 (1996); Boxio, R. et al. Scand. *J. Immunol.* **62**, 140 (2005); Seo, J. et al. *J. Immunol.* **158**, 1895 (1997).

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.

Leukocytes were collected by abdominal lavage 12–14 h after zymosan injection and PMN were isolated by density gradient centrifugation as described above. fMLP (10 nM; Sigma-Aldrich) or WKYMVm (1 nM; AnaSpec) resuspended in Gey's balanced salt solution (Sigma-Aldrich) was added in the lower chamber and 10<sup>6</sup> PMN were incubated in the upper chamber at 37°C in 5% CO<sub>2</sub> for 2 h. After incubation, filters were fixed with methanol and stained with hematoxylin. Cells that migrated to the basal side of the membrane were counted under a microscope and expressed as cells per square millimeter membrane surface-Gyurko, R. et al. J. Immunol. 177, 7250 (2006).

## **Published Citations:**

Gyurko, R. et al. J. Immunol. 177, 7250 (2006).

## Related Products:

Name Cat # Size
WRW4, Formyl Peptide Receptor-Like 1 (FPRL1) Antagonist AS-62638 1 mg
WRWWWW

For Research Use Only