

# PRODUCT INFORMATION



## WKYMVm (trifluoroacetate salt)

Item No. 33589

**Formal Name:** L-tryptophyl-L-lysyl-L-tyrosyl-L-methionyl-L-valyl-D-methioninamide, trifluoroacetate salt

**Synonym:** Trp-Lys-Tyr-Met-Val-D-Met

**MF:** C<sub>41</sub>H<sub>61</sub>N<sub>9</sub>O<sub>7</sub>S<sub>2</sub> • XCF<sub>3</sub>COOH

**FW:** 856.1

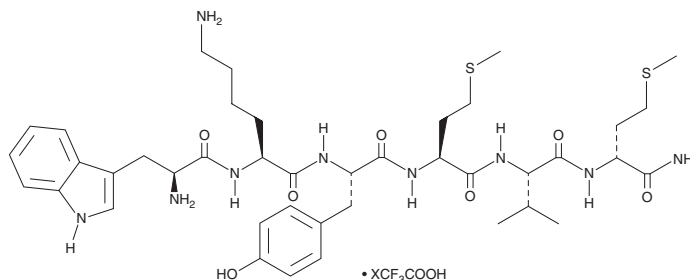
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 219 nm

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

WKYMVm (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the WKYMVm (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. WKYMVm (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of WKYMVm (trifluoroacetate salt) in these solvents is approximately 30 mg/ml.

WKYMVm (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, WKYMVm (trifluoroacetate salt) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. WKYMVm (trifluoroacetate salt) has a solubility of approximately 0.20 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

WKYMVm is a synthetic peptide agonist of formyl peptide receptor 1 (FPR1) and FPR2, which was previously known as formyl peptide receptor-like 1 (FPRL1).<sup>1-3</sup> It induces calcium mobilization in ETFR rat basophilic leukemia cells transfected with FPR1 or FPR2 when used at a concentration of 0.1 nM.<sup>1</sup> WKYMVm stimulates chemotaxis of monocytes, dendritic cells, and natural killer (NK) cells, as well as induces superoxide production in monocytes and neutrophils, *in vitro*.<sup>3</sup> *In vivo*, WKYMVm (8 mg/kg) reverses mucosal destruction, decreases in body weight, and colonic shortening in a mouse model of colitis induced by dextran sulfate (sodium salt) (DSS; Item No. 23250). Topical application of WKYMVm stimulates angiogenesis and accelerates re-epithelialization and granulation tissue formation in a rat model of diabetic cutaneous wounds.<sup>2</sup>

### References

1. Le, Y., Gong, W., Li, B., *et al.* Utilization of two seven-transmembrane, G protein-coupled receptors, formyl peptide receptor-like 1 and formyl peptide receptor, by the synthetic hexapeptide WKYMVm for human phagocyte activation. *J. Immunol.* **163**(12), 6777-6784 (1999).
2. Kwon, Y.W., Heo, S.C., Jang, I.H., *et al.* Stimulation of cutaneous wound healing by an FPR2-specific peptide agonist WKYMVm. *Wound Repair Regen.* **23**(4), 575-582 (2015).
3. Kim, S.D., Kwon, S., Lee, S.K., *et al.* The immune-stimulating peptide WKYMVm has therapeutic effects against ulcerative colitis. *Exp. Mol. Med.* **45**(9), e40 (2013).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/07/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM