

# PRODUCT INFORMATION



## Substance P (7-11) (trifluoroacetate salt)

Item No. 43080

**Formal Name:** L-phenylalanyl-L-phenylalanylglycyl-L-leucyl-L-methioninamide, trifluoroacetate salt

**Synonyms:** Phe-Phe-Gly-Leu-Met, SP (7-11)

**Peptide Sequence:** FFGLM-NH<sub>2</sub>

**MF:** C<sub>31</sub>H<sub>44</sub>N<sub>6</sub>O<sub>5</sub>S • XCF<sub>3</sub>COOH

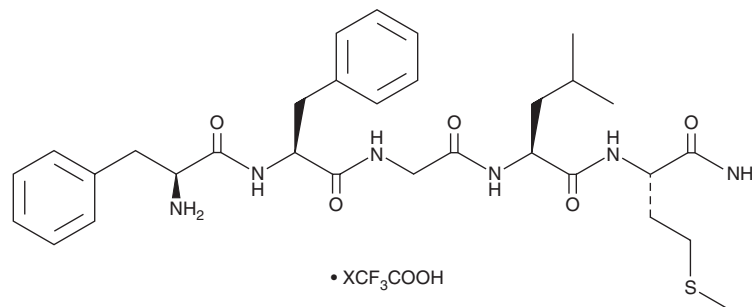
**FW:** 612.8

**Purity:** ≥98%

**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

Substance P (7-11) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the substance P (7-11) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Substance P (7-11) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

### Description

Substance P (7-11) is a peptide fragment of substance P (Item No. 24035).<sup>1</sup> It is formed from substance P via hydrolysis by neutral endopeptidase. Substance P (7-11) (>1 μM) increases collagenase and prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14041) production in isolated bovine articular chondrocytes.<sup>2</sup> It increases intracellular calcium levels in isolated bovine chondrocytes when used at concentrations ranging from 0.01 to 10 μM.<sup>1</sup>

### References

1. Halliday, D.A., McNeil, J.D., and Scicchitano, R. A metabolite of substance P, SP<sub>7-11</sub> is involved in the pathogenesis of inflammatory joint disease. *Med. Hypotheses* **40(4)**, 227-231 (1993).
2. Halliday, D.A., McNeil, J.D., Betts, W.H., et al. The substance P fragment SP-(7-11) increases prostaglandin E<sub>2</sub>, intracellular Ca<sup>2+</sup> and collagenase production in bovine articular chondrocytes. *Biochem. J.* **292(1)**, 57-62 (1993).

#### 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, 12/03/2025

#### 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