

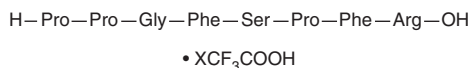
# PRODUCT INFORMATION



## Bradykinin (2-9) (trifluoroacetate salt)

Item No. 39723

<b>Synonyms:</b>	BK (2-9), (Des-Arg <sup>1</sup> )-BK, (Des-Arg <sup>1</sup> )-Bradykinin, H-Pro-Pro-Gly-Phe-Ser-Pro-Phe-Arg-OH
<b>Peptide Sequence:</b>	PPGFSPFR-OH
<b>MF:</b>	C <sub>44</sub> H <sub>61</sub> N <sub>11</sub> O <sub>10</sub> • XCF <sub>3</sub> COOH
<b>FW:</b>	904.0
<b>Purity:</b>	≥98%
<b>Supplied as:</b>	A solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥4 years



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

### Laboratory Procedures

Bradykinin (2-9) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the bradykinin (2-9) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Bradykinin (2-9) is a metabolite of the endogenous vasodilator peptide bradykinin (Item No. 37408).<sup>1</sup> It is formed from bradykinin via N-terminal cleavage by carboxypeptidase.

### Reference

1. Gangnus, T. and Burckhardt, B.B. Targeted LC-MS/MS platform for the comprehensive determination of peptides in the kallikrein-kinin system. *Anal. Bioanal. Chem.* **413(11)**, 2971-2984 (2023).

#### 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, 02/22/2024

#### 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