

PRODUCT INFORMATION

Lys-(Des-Arg⁹, Leu⁸)-Bradykinin (trifluoroacetate salt)

Item No. 35462

Formal Name:	9-L-leucine-1-9-kallidin, trifluoroacetate salt		
Synonyms:	[Des-Arg ¹⁰ ,Leu ⁹]-Kallidin, [Des-Arg ¹⁰ ,Leu ⁹]-KD, Lys[Leu ⁸]Des-Arg ⁹ -BK, Lys[Leu ⁸]Des-Arg ⁹ -Bradykinin		
MF:	C ₄₇ H ₇₅ N ₁₃ O ₁₁ • XCF ₃ COOH	H-Lys-Arg-Pro-Pro-Gly-Phe-Ser-Pro-Leu-OH	• XCF ₃ COOH
FW:	998.2		
Purity:	≥95%		
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

Lys-(Des-Arg⁹, Leu⁸)-Bradykinin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Lys-(Des-Arg⁹, Leu⁸)-bradykinin (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Lys-(Des-Arg⁹, Leu⁸)-Bradykinin (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Lys-(Des-Arg⁹, Leu⁸)-bradykinin (trifluoroacetate salt) in these solvents is approximately 3 and 1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Lys-(Des-Arg⁹, Leu⁸)-bradykinin (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of Lys-(Des-Arg⁹, Leu⁸)-bradykinin (trifluoroacetate salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Lys-(Des-Arg⁹, Leu⁸)-Bradykinin is a bradykinin B₁ receptor antagonist (K_i = 0.43 nM for the rabbit aorta receptor).¹ It is selective for the bradykinin B₁ over the B₂ receptor (K_i = >10 μM for the rabbit ileum receptor).² Lys-(Des-Arg⁹, Leu⁸)-Bradykinin inhibits contractions induced by the bradykinin B₁ receptor agonist Des-Arg⁹-bradykinin (Item No. 31136) in isolated rabbit aortic rings (pA₂ = 8.54).³ Intraperitoneal administration of Lys-(Des-Arg⁹, Leu⁸)-bradykinin (5 mg/kg per day for three days) reduces increases in paw edema in a rat model of arthritis induced by peptidoglycan-polysaccharide (PGPS).⁴

References

- MacNeil, T., Bierilo, K.K., Menke, J.G., et al. Cloning and pharmacological characterization of a rabbit bradykinin B₁ receptor. *Biochimica et Biophysica Acta* **1264**(2), 223-228 (1995).
- Meini, S., Cucchi, P., Catalani, C., et al. Radioligand binding characterization of the bradykinin B₂ receptor in the rabbit and pig ileal smooth muscle. *Eur. J. Pharmacol.* **635**(1-3), 34-39 (2010).
- Drapeau, G., Audet, R., Levesque, L., et al. Development and *in vivo* evaluation of metabolically resistant antagonists of B₁ receptors for kinins. *J. Pharmacol. Exp. Ther.* **266**(1), 192-199 (1993).
- Blais, C., Jr., Couture, R., Drapeau, G., et al. Involvement of endogenous kinins in the pathogenesis of peptidoglycan-induced arthritis in the Lewis rat. *Arthritis Rheum.* **40**(7), 1327-1333 (1997).

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, 11/14/2022

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