PRODUCT INFORMATION



Spinorphin

Item No 29914

CAS Registry No.:	137201-62-8	N-H
Formal Name:	L-leucyl-L-valyl-L-valyl-L-tyrosyl-L-	
	prolyl-L-tryptophyl-L-threonine	N N OH
Synonyms:	Leu-Val-Val-Tyr-Pro-Trp-Thr,	
	LVVYPWT	
MF:	$C_{45}H_{44}N_{0}O_{10}$	
FW:	877.0	
Purity:	≥95%	
UV/Vis.:	λ: 222 nm	
Supplied as:	A crystalline solid	
Storage:	-20°C	о́н
Stability:	≥4 years	
-	•	

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

Laboratory Procedures

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

Description

Spinorphin is a heptapeptide inhibitor of the enkephalin-degrading enzymes aminopeptidase, dipeptidyl aminopeptidase, angiotensin-converting enzyme (ACE), and enkephalinase (IC₅₀s = 3.3, 1.4, 2.4, and 10 µg/ml, respectively, for monkey brain enzymes).¹ It is selective for these enzymes over human serum aminopeptidase A (IC₅₀ = >100 μ g/ml), as well as porcine kidney aminopeptidase B, aminopeptidase M, dipeptidyl peptidase 1 (DPP-1), DPP-2, DPP-3, and DPP-4 (IC₅₀s = >55 μ g/ml for all). Spinorphin inhibits chemotaxis, production of reactive oxygen species (ROS), and exocytosis of glucuronidase and collagenase in polymorphonuclear neutrophils (PMNs). It potentiates enkephalin-induced action potentials in rat hippocampal slices. Spinorphin (5 µg/animal) enhances the antinociceptive effects of Leu-enkephalin (Item No. 23283) in mice.²

References

- 1. Yamamoto, Y., Ono, H., Ueda, A., et al. Spinorphin as an endogenous inhibitor of enkephalin-degrading enzymes: Roles in pain and inflammation. Curr. Protein Pept. Sci. 3(6), 587-599 (2002).
- 2. Honda, M., Okutsu, H., Matsuura, T., et al. Spinorphin, an endogenous inhibitor of enkephalin-degrading enzymes, potentiates leu-enkephalin-induced anti-allodynic and antinociceptive effects in mice. Jpn. J. Pharmacol. 87(4), 261-267 (2001).

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

SAFETY DATA

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

uyer 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, 12/13/2022

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