

PRODUCT INFORMATION



Lysipressin (acetate)

Item No. 37430

CAS Registry No.: 83968-49-4
Formal Name: 8-L-lysine-vasopressin, monoacetate
Synonyms: CYFQNCPKG-NH₂, Lypressin, Lys⁸-vasopressin, Lysine⁸-vasopressin, Lysine Vasopressin

MF: C₄₆H₆₅N₁₃O₁₂S₂ • C₂H₄O₂

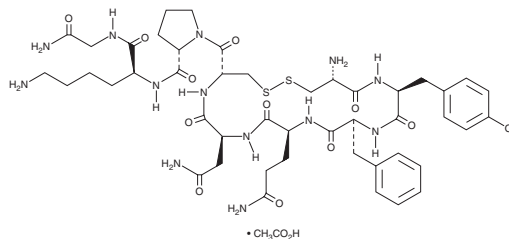
FW: 1,116.3

Purity: ≥98%

Supplied as: A solid

Storage: -80°C

Stability: ≥4 years



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

Laboratory Procedures

Lysipressin (acetate) is supplied as a solid. A stock solution may be made by dissolving the lysipressin (acetate) in the solvent of choice, which should be purged with an inert gas. Lysipressin (acetate) is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

Lysipressin is a peptide hormone found in pigs and marsupials that corresponds to amino acids 20-28 of the vasoconstrictor peptide vasopressin with lysine in place of arginine at position 27.^{1,2} It is an arginine vasopressin receptor 1A (V_{1A}) receptor agonist that binds to rat and human V_{1A} receptors (K_s = 1.2 and 10.1 nM, respectively) and increases intracellular calcium levels in cultured rat and human aortic cells with half-maximal activation (K_{act}) values of 12 and 20.8 nM, respectively.³ Lysipressin induces vasoconstriction in isolated mesenteric arteries from normotensive and spontaneously hypertensive rats.⁴ It also has antidiuretic effects in non-anesthetized dwarf pigs, decreasing urine volume and increasing urinary solutes without affecting the glomerular filtration rate.¹

References

1. Munsick, R.A., Sawyer, W.H., and Van Dyke, H.B. The antidiuretic potency of arginine and lysine vasopressins in the pig with observations on porcine renal function. *Endocrinology* **63**(5), 688-693 (1958).
2. Chauvet, J., Hurpet, D., Michel, G., et al. Two multigene families for marsupial neurohypophysial hormones? Identification of oxytocin, mesotocin, lysipressin and arginine vasopressin in the North American opossum (*Didelphis virginiana*). *Biochem. Biophys. Res. Commun.* **123**(1), 306-311 (1984).
3. Colson, P.H., Virsolvy, A., Gaudard, P., et al. Terlipressin, a vasoactive prodrug recommended in hepatorenal syndrome, is an agonist of human V1, V2 and V1B receptors: Implications for its safety profile. *Pharmacol. Res.* **113**(Pt A), 257-264 (2016).
4. Schlegel, P.A., Monney, M., and Brunner, H.R. Isolated perfused mesenteric arteries of hypertensive and normotensive rats; response to norepinephrine, lysine vasopressin and angiotensin II. *Clin. Exp. Hypertens. A.* **7**(11), 1583-1596 (1985).

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, 04/10/2026

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