

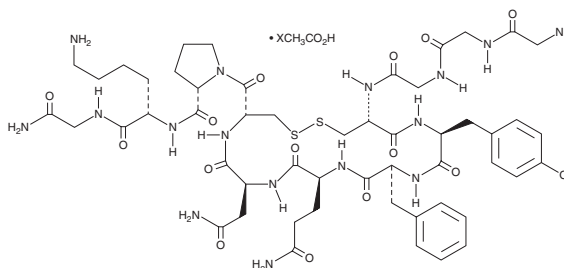
PRODUCT INFORMATION



Terlipressin (acetate)

Item No. 27027

CAS Registry No.: 914453-96-6
Formal Name: N-(glycylglycylglycyl)-8-L-lysine-vasopressin, acetate
Synonyms: Glycylpressin, Remestyp
MF: $C_{52}H_{74}N_{16}O_{15}S_2 \cdot XC_2H_4O_2$
FW: 1,227.4
Purity: $\geq 95\%$
Supplied as: A crystalline 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

Terlipressin (acetate) is supplied as a crystalline solid. A stock solution may be made by dissolving the terlipressin (acetate) in the solvent of choice, which should be purged with an inert gas. Terlipressin (acetate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of terlipressin (acetate) in these solvents is approximately 25 and 20 mg/ml, respectively. Terlipressin (acetate) is also slightly soluble in ethanol.

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 terlipressin (acetate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of terlipressin (acetate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Terlipressin is a prodrug form of the vasopressin peptide lysipressin and a partial agonist of the vasopressin V_{1A} receptor ($K_i = 0.85 \mu\text{M}$).¹ It is also an agonist of vasopressin V_{1B} and V_2 receptors ($K_s = 1.11$ and $1.58 \mu\text{M}$, respectively). It increases mean arterial pressure (MAP) and decreases mortality in a rat model of uncontrolled hemorrhagic shock when administered at a dose of $15 \mu\text{g/kg}$.² Terlipressin ($2.6 \mu\text{g/kg}$ per hour), in combination with norepinephrine, improves vascular reactivity and increases survival time in a rat model of cecal ligation and puncture-induced septic shock and a rabbit model of LPS-induced septic shock.³ It also increases MAP in a rat model of liver cirrhosis with portal hypertension, arterial hypotension, high cardiac output, and low systemic vascular resistance.⁴

References

1. Colson, P.H., Virsolvy, A., Gaudard, P., *et al.* Terlipressin, a vasoactive prodrug recommended in hepatorenal syndrome, is an agonist of human V_1 , V_2 and V_{1B} receptors: Implications for its safety profile. *Pharmacol. Res.* **113(Pt A)**, 257-264 (2016).
2. Lee, C.C., Lee, M.T., Chang, S.S., *et al.* A comparison of vasopressin, terlipressin, and lactated ringers for resuscitation of uncontrolled hemorrhagic shock in an animal model. *PLoS One* **9(4)**, e95821 (2014).
3. Xiao, X., Zhu, Y., Zhen, D., *et al.* Beneficial and side effects of arginine vasopressin and terlipressin for septic shock. *J. Surg. Res.* **195(2)**, 568-579 (2015).
4. Fernández-Varo, G., Oró, D., Cable, E.E., *et al.* Vasopressin $1a$ receptor partial agonism increases sodium excretion and reduces portal hypertension and ascites in cirrhotic rats. *Hepatology* **63(1)**, 207-216 (2016).

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, 11/07/2022

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