PRODUCT INFORMATION



Angiotensin II (human) (acetate)

Item No. 17150

		NH		
CAS Registry No.:	68521-88-0	N	○ 0 ^H	H /
Formal Name:	5-L-isoleucine-angiotensin II, acetate	н		N
MF:	$C_{50}H_{71}N_{13}O_{12} \bullet XC_{2}H_{4}O_{2}$			N O OH
FW:	1,046.2	0 ^H 0	H O	
Purity:	≥90%	HO	N N N N	
Supplied as:	A crystalline solid	Ü A _{H₂} Ü ¦	L U L	
Storage:	-20°C	/ \		~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~~
Stability:	≥4 years	•	XCH ₃ CO ₂ H	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis				

нм

Laboratory Procedures

Angiotensin II (human) (acetate) is supplied as a crystalline solid. Aqueous solutions of angiotensin II (human) (acetate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of angiotensin II (human) (acetate) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Angiotensin II is a peptide hormone known best as a vasoconstrictor with central roles in chronic hypertension, heart failure, and stroke.¹ It is an octapeptide typically generated by the removal of two residues from angiotensin I by angiotensin-converting enzyme (ACE).² Angiotensin II is a ligand for at least two distinct receptors, AT_1 and AT_2 , each evoking distinct signaling pathways and physiological responses.^{1,3} The development of antagonists for specific angiotensin II receptor subtypes represents a valuable alternative to ACE inhibitors.⁴

References

- 1. Burnier, M. Angiotensin II type 1 receptor blockers. Circulation 103(6), 904-912 (2001).
- 2. Redelinghuys, P., Nchinda, A.T., and Sturrock, E.D. Development of domain-selective angiotensin I-converting enzyme inhibitors. Ann. N. Y. Acad. Sci. 1056, 160-175 (2005).
- 3. Bosnyak, S., Jones, E.S., Christopoulos, A., et al. Relative affinity of angiotensin peptides and novel ligands at AT₁ and AT₂ receptors. Clin. Sci. (Lond) **121(7)**, 297-303 (2011).
- Wexler, R.R., Greenlee, W.J., Irvin, J.D., et al. Nonpeptide angiotensin II receptor antagonists: The next 4. generation in antihypertensive therapy. J. Med. Chem. 39(3), 625-656 (2014).

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

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, 10/25/2022

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