

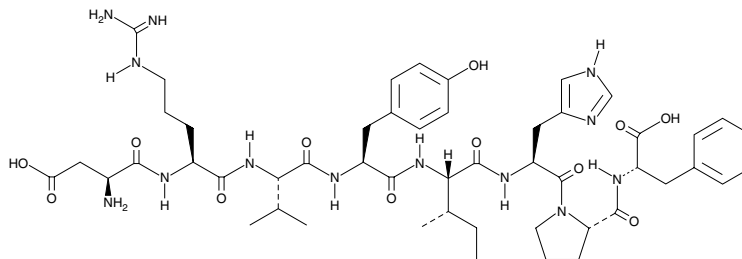
# Product Information



## Angiotensin II (human)

Item No. 17150

**CAS Registry No.:** 4474-91-3  
**Formal Name:** 5-L-isoleucine-angiotensin II  
**MF:** C<sub>50</sub>H<sub>71</sub>N<sub>13</sub>O<sub>12</sub>  
**FW:** 1,046.2  
**Purity:** ≥90%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that angiotensin II (human) be stored as supplied at -20°C. It should be stable for at least two years.

Angiotensin II (human) is supplied as a crystalline solid. Aqueous solutions of angiotensin II (human) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of angiotensin II (human) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

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

### References

1. Burnier, M. Angiotensin II type 1 receptor blockers. *Circulation* **103**, 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 AT<sub>1</sub> and AT<sub>2</sub> receptors. *Clin. Sci. (Lond)* **121**(7), 297-303 (2011).
4. Wexler, R.R., Greenlee, W.J., Irvin, J.D., *et al.* Nonpeptide angiotensin II receptor antagonists: The next generation in antihypertensive therapy. *J. Med. Chem.* **39**(3), 625-656 (2014).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/17150](http://www.caymanchem.com/catalog/17150)

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
[www.caymanchem.com](http://www.caymanchem.com)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 02/13/2015