

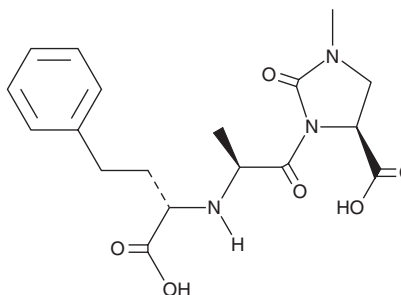
# PRODUCT INFORMATION



## Imidaprilat

Item No. 38046

**CAS Registry No.:** 89371-44-8  
**Formal Name:** 3-[(2S)-2-[[[(1S)-1-carboxy-3-phenylpropyl]amino]-1-oxopropyl]-1-methyl-2-oxo-4S-imidazolidinecarboxylic acid  
**Synonym:** 6366A  
**MF:** C<sub>18</sub>H<sub>23</sub>N<sub>3</sub>O<sub>6</sub>  
**FW:** 377.4  
**Purity:** ≥95%  
**Supplied as:** A 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

Imidaprilat is supplied as a solid. A stock solution may be made by dissolving the imidaprilat in the solvent of choice, which should be purged with an inert gas. Imidaprilat is soluble in the organic solvent DMSO.

### Description

Imidaprilat is an inhibitor of angiotensin-converting enzyme (ACE; IC<sub>50</sub> = 2.6 nM) and an active metabolite of the prodrug imidapril (Item No. 18854).<sup>1</sup> It inhibits bradykinin-induced contractions in isolated guinea pig ileum (IC<sub>50</sub> = 1.7 nM). Imidaprilat (50 nM) reduces adhesion induced by chemokine (C-C-) motif ligand 2 (CCL2) of THP-1 monocytes to TNF-α-activated human umbilical vein endothelial cells (HUVECs).<sup>2</sup> *In vivo*, imidaprilat (1 mg/kg) inhibits the angiotensin I-induced pressor response in unanesthetized normotensive rats.<sup>1</sup>

### References

1. Kubo, M., Kato, J., Ochiai, T., *et al.* Pharmacological studies on (4S)-1-methyl-3-[(2S)-2-[N-((1S)-1-ethoxycarbonyl-3-phenylpropyl)amino] propionyl]-2-oxo-imidazolidine-4-carboxylic acid hydrochloride (TA-6366), a new ACE inhibitor: I. ACE inhibitory and anti-hypertensive activities. *Jpn. J. Pharmacol.* **53**(2), 201-210 (1990).
2. Kojima, C., Kawakami, A., Takei, T., *et al.* Angiotensin-converting enzyme inhibitor attenuates monocyte adhesion to vascular endothelium through modulation of intracellular zinc. *J. Pharmacol. Exp. Ther.* **323**(3), 855-860 (2007).

#### 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, 02/12/2024

#### 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