

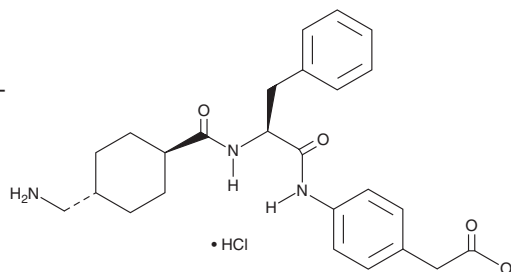
# PRODUCT INFORMATION



## PKSI-527

Item No. 23974

**CAS Registry No.:** 128837-71-8  
**Formal Name:** 4-[[[(2S)-2-[[[trans-4-(aminomethyl)cyclohexyl]carbonyl]amino]-1-oxo-3-phenylpropyl]amino]-benzeneacetic acid, monohydrochloride  
**MF:** C<sub>25</sub>H<sub>31</sub>N<sub>3</sub>O<sub>4</sub> • HCl  
**FW:** 474.0  
**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

PKSI-527 is supplied as a solid. A stock solution may be made by dissolving the PKSI-527 in water. The solubility of PKSI-527 in water is approximately 10 mM. We do not recommend storing the aqueous solution for more than one day.

### Description

PKSI-527 is an inhibitor of plasma kallikrein ( $K_i = 0.81 \mu\text{M}$ ).<sup>1</sup> It is selective for plasma kallikrein over glandular kallikrein, plasmin, thrombin, urokinase, and Factor Xa ( $K_i$ s = >500, 390, >500, 200, and >500  $\mu\text{M}$ , respectively). PKSI-527 reduces bradykinin generation induced by kaolin and  $\lambda$ -carrageenan *ex vivo* in human plasma. It also prolongs partial thromboplastin and euglobulin clot lysis times. *In vivo*, PKSI-527 (300 mg/kg per day) reduces hyperplasia, pannus formation, and infiltration of inflammatory cells in the tarsal joint of mice with collagen-induced arthritis.<sup>2</sup>

### References

1. Wanaka, K., Okamoto, S., Bohgaki, M., *et al.* Effect of a highly selective plasma-kallikrein synthetic inhibitor on contact activation relating to kinin generation, coagulation and fibrinolysis. *Thromb. Res.* **57(6)**, 889-895 (1990).
2. Fujimori, Y., Nakamura, T., Shimizu, K., *et al.* Effects of a highly selective plasma kallikrein inhibitor on collagen-induced arthritis in mice. *Agents Actions* **39(1-2)**, 42-48 (1993).

#### 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, 12/14/2022

#### 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