

PRODUCT INFORMATION

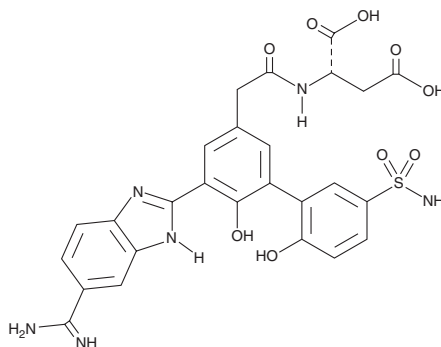


PCI 27483

Item No. 21334

CAS Registry No.: 871266-63-6
Formal Name: N-[2-[5-[6-(aminoiminomethyl)-1H-benzimidazol-2-yl]-5'-(aminosulfonyl)-2',6'-dihydroxy[1,1'-biphenyl]-3-yl]acetyl]-L-aspartic acid

MF: C₂₆H₂₄N₆O₉S
FW: 596.6
Purity: ≥95%
UV/Vis.: λ_{max}: 230, 307, 340 nm
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

PCI 27483 is supplied as a crystalline solid. A stock solution may be made by dissolving the PCI 27483 in the solvent of choice. PCI 27483 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of PCI 27483 in these solvents is approximately 10 and 1 mg/ml, respectively.

PCI 27483 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PCI 27483 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PCI 27483 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PCI 27483 is a small molecule inhibitor of activated coagulation factor VIIa with antithrombotic effects in a baboon model of arterial thrombosis.¹ It also inhibits the growth of BxPC3 xenografts by 42 and 85% at doses of 60 and 90 mg/kg, respectively.²

References

1. Gómez-Outes, A., Suárez-Gea, L., Lecumberri, R., *et al.* New parenteral anticoagulants in development. *Ther. Adv. Cardiovasc. Disc.* **5**(1), 33-59 (2011).
2. Prescott, J., Thiemann, P., and Loury, D. PCI-27483, a small molecule inhibitor of Factor VIIa, inhibits growth of BxPC3 pancreatic adenocarcinoma xenograft tumors. *AACR Annual Meeting* (2008).

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, 10/24/2022

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