

PRODUCT INFORMATION

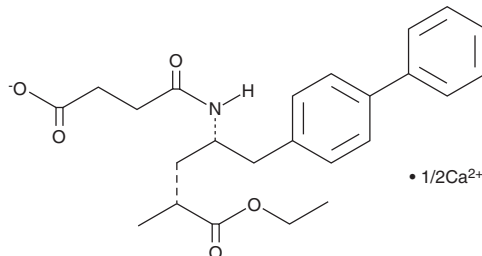


AHU377 (calcium salt)

Item No. 21473

CAS Registry No.: 1369773-39-6
Formal Name: (αR,γS)-γ-[(3-carboxy-1-oxopropyl)amino]-α-methyl-[1,1'-biphenyl]-4-pentanoic acid, 4-ethyl ester, hemicalcium salt

Synonym: Sacubitril
MF: C₂₄H₂₈NO₅ • 1/2Ca
FW: 430.5
Purity: ≥98%
UV/Vis.: λ_{max}: 254 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

AHU377 (calcium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the AHU377 (calcium salt) in the solvent of choice, which should be purged with an inert gas. AHU377 (calcium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AHU377 (calcium salt) in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

AHU377 (calcium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AHU377 (calcium salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AHU377 (calcium salt) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AHU377 is an ethyl ester prodrug from of the neprilysin inhibitor LBQ657 (Item No. 19829).¹ It increases plasma levels of atrial natriuretic factor (ANF) in conscious rats and anesthetized dogs when administered at doses of 10 and 30 mg/kg, respectively. AHU377 (30 mg/kg) increases ANF-induced urinary sodium excretion in anesthetized rats. Formulations containing AHU377 in combination with valsartan have been used for the treatment of mild to moderate hypertension and chronic heart failure.

Reference

1. Ksander, G.M., Ghai, R.D., deJesus, R., *et al.* Dicarboxylic acid dipeptide neutral endopeptidase inhibitors. *J. Med. Chem.* **38**(10), 1689-1700 (1995).

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, 12/13/2022

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