

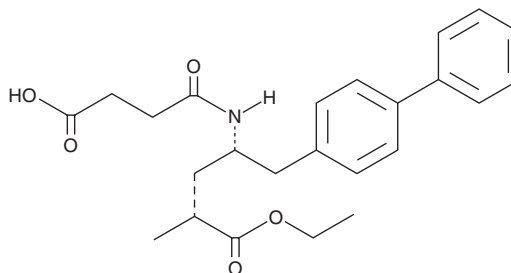
PRODUCT INFORMATION



AHU377

Item No. 36260

CAS Registry No.: 149709-62-6
Formal Name: (αR,γS)-γ-[(3-carboxy-1-oxopropyl)amino]-α-methyl-[1,1'-biphenyl]-4-pentanoic acid, α-ethyl ester
Synonym: Sacubitril
MF: C₂₄H₂₉NO₅
FW: 411.5
Purity: ≥98%
UV/Vis.: λ_{max}: 252 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

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 in 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, 12/12/2022

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