

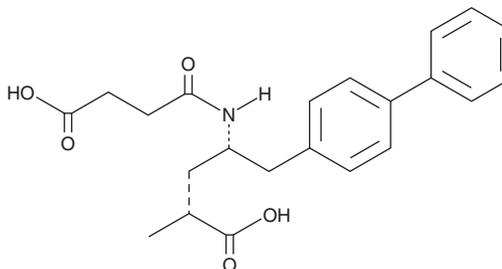
PRODUCT INFORMATION



LBQ657

Item No. 19829

CAS Registry No.: 149709-44-4
Formal Name: γ S-[(3-carboxy-1-oxopropyl) amino]- α R-methyl-[1,1'-biphenyl]-4-pentanoic acid
Synonym: Sacubitrilat
MF: C₂₂H₂₅NO₅
FW: 383.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 254 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

LBQ657 is supplied as a crystalline solid. A stock solution may be made by dissolving the LBQ657 in the solvent of choice. LBQ657 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of LBQ657 in these solvents is approximately 30 mg/ml.

LBQ657 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LBQ657 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. LBQ657 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LBQ657 is an inhibitor of neprilysin ($IC_{50} = 5 \text{ nM}$).¹ It is selective for neprilysin over angiotensin converting enzyme (ACE) when used at a concentration of 10 μM . LBQ657 inhibits neprilysin cleavage of enkephalin, glutaryl-Ala-Ala-Phe- β -naphthylamide, and atrial natriuretic factor (IC_{50} s = 5, 0.74, and 5.3 nM, respectively).

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, 03/27/2025

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