

# PRODUCT INFORMATION

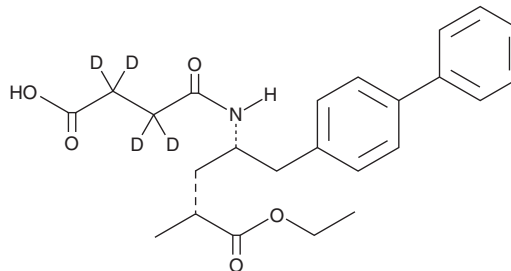


AHU377-d<sub>4</sub>  
Item No. 25264

**CAS Registry No.:** 1884269-07-1  
**Formal Name:** (αR,γS)-γ-[(3-carboxy-1-oxopropyl-2,2,3,3-d<sub>4</sub>)amino]-α-methyl-[1,1'-biphenyl]-4-pentanoic acid, α-ethyl ester

**Synonym:** Sacubitril-d<sub>4</sub>  
**MF:** C<sub>24</sub>H<sub>25</sub>D<sub>4</sub>NO<sub>5</sub>  
**FW:** 415.5  
**Chemical Purity:** ≥98% (AHU377)

**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>4</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A film  
**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-d<sub>4</sub> is intended for use as an internal standard for the quantification of AHU377 (Item Nos. 36260 | 21473) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

AHU377-d<sub>4</sub> is supplied as a film. A stock solution may be made by dissolving the AHU377-d<sub>4</sub> in the solvent of choice, which should be purged with an inert gas. AHU377-d<sub>4</sub> is slightly soluble in chloroform, DMSO, and methanol.

## Description

AHU377 is an ethyl ester prodrug form of the neprilysin inhibitor LBQ657 (Item No. 19829).<sup>1</sup> 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