

# PRODUCT INFORMATION



## LHW090-A7

Item No. 18410

**Formal Name:** (αS)-3'-chloro-α-[[[(1S)-2-ethoxy-1-methyl-2-oxoethyl]amino]-[1,1'-biphenyl]-4-propanoic acid, monosodium salt

**MF:** C<sub>20</sub>H<sub>21</sub>ClNO<sub>4</sub> • Na

**FW:** 397.8

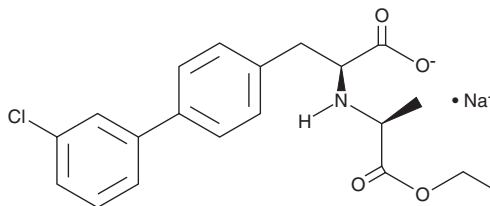
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 257 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

LHW090-A7 is supplied as a crystalline solid. A stock solution may be made by dissolving the LHW090-A7 in the solvent of choice, which should be purged with an inert gas. LHW090-A7 is soluble in organic solvents such as ethanol and DMSO. The solubility of LHW090-A7 in these solvents is approximately 1 and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of LHW090-A7 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of LHW090-A7 in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

LHW090-A7 is a compound developed as an inhibitor of neutral endopeptidase (NEP), an enzyme responsible for the metabolic inactivation of enkephalins.<sup>1</sup> Inhibitors of NEP have been shown to increase plasma levels of endogenous atrial natriuretic peptides and, as such, are expected to induce natriuretic and diuretic effects.

### Reference

1. Iwaki, Y., Kawanami, T., Ksander, G.M., *et al.* Substituted carbamoylmethylamino acetic acid derivatives as novel NEP inhibitors. WO 2011/061271 A1 (2011), PCT/EP2010/067781, Novartis AG.

#### 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