

# PRODUCT INFORMATION



## Encaleret

Item No. 45051

**CAS Registry No.:** 787583-71-5  
**Formal Name:** 2'-[(1R)-1-[(2R)-3-[[2-(4-chloro-3-fluorophenyl)-1,1-dimethylethyl]amino]-2-hydroxypropoxy]ethyl]-3-methyl-[1,1'-biphenyl]-4-carboxylic acid

**Synonyms:** CLTX-305, JTT-305, MK-5442

**MF:** C<sub>29</sub>H<sub>33</sub>ClFNO<sub>4</sub>

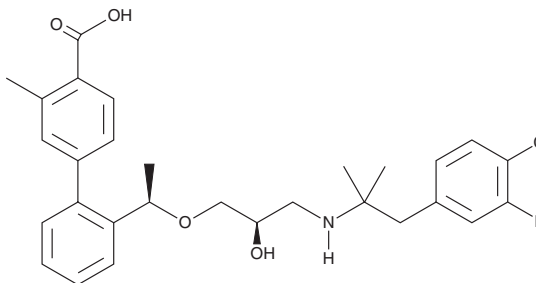
**FW:** 514.0

**Purity:** ≥95%

**Supplied as:** A 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

Encaleret is supplied as a solid. A stock solution may be made by dissolving the encaleret in the solvent of choice, which should be purged with an inert gas. Encaleret is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

### Description

Encaleret is a calcium-sensing receptor (CaSR) antagonist (IC<sub>50</sub> = 0.012 μM in a reporter assay using PC12h cells).<sup>1</sup> Oral administration of encaleret (20 μg/g) increases plasma levels of parathyroid hormone, prevents renal calcification, and increases bone mineral density in a model of autosomal dominant hypocalcemia (ADH) using CaSR knock-in mice.<sup>2</sup>

### References

1. Shinagawa, Y., Inoue, T., Katsushima, T., *et al.* Discovery of a potent and short-acting oral calcilytic with a pulsatile secretion of parathyroid hormone. *ACS Med. Chem. Lett.* **2(3)**, 238-242 (2010).
2. Dong, B., Endo, I., Ohnishi, Y., *et al.* Calcilytic ameliorates abnormalities of mutant calcium-sensing receptor (CaSR) knock-in mice mimicking autosomal dominant hypocalcemia (ADH). *J. Bone Miner. Res.* **30(11)**, 1980-1993 (2015).

#### 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, 02/13/2026

#### 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](http://WWW.CAYMANCHEM.COM)