

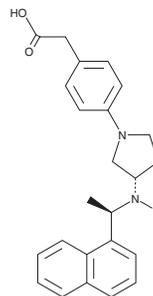
PRODUCT INFORMATION



Evocalcet

Item No. 33346

CAS Registry No.: 870964-67-3
Formal Name: 4-[(3S)-3-[[[(1R)-1-(1-naphthalenyl)ethyl]amino]-1-pyrrolidinyl]-benzeneacetic acid
Synonyms: KHK7580, MT-4580
MF: C₂₄H₂₆N₂O₂
FW: 374.5
Purity: ≥98%
UV/Vis.: λ_{max}: 224 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

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

Evocalcet is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, evocalcet should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Evocalcet has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Evocalcet is an agonist of the calcium-sensing receptor (CaSR).¹ It induces intracellular calcium release in HEK293 cells expressing human CaSR (EC₅₀ = 92.7 nM). Evocalcet reduces serum levels of parathyroid hormone (PTH) and calcium when administered at doses of 0.3 and 0.1 mg/kg, respectively, in non-nephrectomized and 5/6 nephrectomized rats. It also reduces serum levels of PTH without inducing emesis in marmosets when administered at doses of 0.3 and 0.5 mg/kg.

Reference

1. Kawata, T., Tokunaga, S., Murai, M., *et al.* A novel calcimimetic agent, evocalcet (MT4580/KHK7580), suppresses the parathyroid cell function with little effect on the gastrointestinal tract or CYP isozymes *in vivo* and *in vitro*. *PLoS One* **13**(4), e0195316 (2018).

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, 11/14/2022

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