

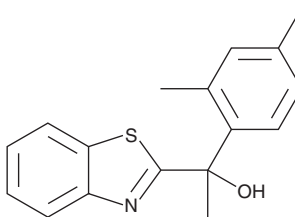
PRODUCT INFORMATION



AC-265347

Item No. 30937

CAS Registry No.: 1253901-26-6
Formal Name: α -(2,4-dimethylphenyl)- α -methyl-2-benzothiazolemethanol
MF: C₁₇H₁₇NOS
FW: 283.4
Purity: \geq 98%
Supplied as: A 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

AC-265347 is supplied as a solid. A stock solution may be made by dissolving the AC-265347 in the solvent of choice, which should be purged with an inert gas. AC-265347 is soluble in organic solvents such as ethanol and DMSO. The solubility of AC-265347 in these solvents is approximately 100 mM.

Description

AC-265347 is a calcimimetic and positive allosteric modulator of the calcium-sensing receptor (CaSR; EC₅₀ = 7.94 nM in a cell-based phosphatidylinositol hydrolysis assay in the presence of calcium).^{1,2} It inhibits the proliferation of NIH3T3 cells expressing human CaSR (EC₅₀ = 25.12 nM).¹ AC-265347 reduces plasma parathyroid hormone levels in rats (ED₅₀ = 0.01 mg/kg).

References

1. Ma, J.N., Owens, M., Gustafsson, M., *et al.* Characterization of highly efficacious allosteric agonists of the human calcium-sensing receptor. *J. Pharmacol. Exp. Ther.* **337**(1), 275-284 (2011).
2. Leach, K., Gregory, K.J., Kufareva, I., *et al.* Towards a structural understanding of allosteric drugs at the human calcium-sensing receptor. *Cell Res.* **26**(5), 574-592 (2016).

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, 11/04/2022

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