

PRODUCT INFORMATION

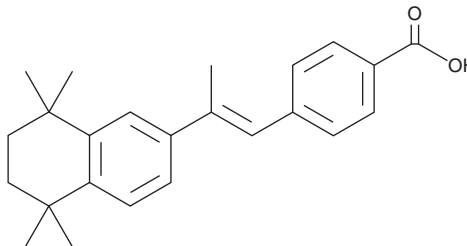


TTNPB

Item No. 16144

CAS Registry No.: 71441-28-6
Formal Name: 4-[(1E)-2-(5,6,7,8-tetrahydro-5,5,8,8-tetramethyl-2-naphthalenyl)-1-propen-1-yl]-benzoic acid
Synonyms: AGN 191183, Arotinoid Acid, Ro 13-7410

MF: C₂₄H₂₈O₂
FW: 348.5
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 303 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

TTNPB is supplied as a crystalline solid. A stock solution may be made by dissolving the TTNPB in the solvent of choice. TTNPB is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of TTNPB in these solvents is approximately 0.1, 2, and 5 mg/ml, respectively.

TTNPB is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TTNPB should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TTNPB has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TTNPB is an analog of retinoic acid (Item No. 11017) that potently and selectively activates retinoic acid receptors (EC₅₀s = 21, 4, and 2.4 nM for RARα, RARβ, and RARγ, respectively).^{1,2} It does not act on retinoid X receptors and weakly agonizes farnesoid X receptor (EC₅₀ > 1 μM).^{2,3} TTNPB is used to study RAR action in diverse processes, including epidermal cell proliferation, embryogenesis, and stem cell differentiation.⁴⁻⁷

References

1. Beard, R.L., Chandraratna, R.A., Colon, D.F., *et al. J. Med. Chem.* **38(15)**, 2820-2829 (1995).
2. Wong, M.F., Repa, J.J., Clagett-Dame, M., *et al. Bioorg. Med. Chem. Lett.* **7(17)**, 2313-2318 (1997).
3. Maloney, P.R., Parks, D.J., Haffner, C.D., *et al. J. Med. Chem.* **43(16)**, 2971-2974 (2000).
4. Minucci, S., Saint-Jeannet, J.-P., Toyama, R., *et al. Proc. Natl. Acad. Sci. USA* **93(5)**, 1803-1807 (1996).
5. Thacher, S.M., Standeven, A.M., Athanikar, J., *et al. J. Pharmacol. Exp. Ther.* **282(2)**, 528-534 (1997).
6. Araoka, T., Mae, S., Kurose, Y., *et al. PLoS One* **9(1)**, e84881 (2014).
7. Hou, P., Li, Y., Zhang, X., *et al. Science* **341(6146)**, 651-654 (2013).

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, 06/25/2020

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