

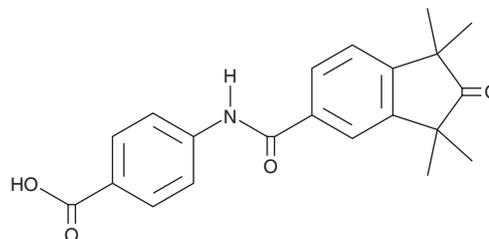
PRODUCT INFORMATION



BMS 753

Item No. 29493

CAS Registry No.: 215307-86-1
Formal Name: 4-[[[(2,3-dihydro-1,1,3,3-tetramethyl-2-oxo-1H-inden-5-yl)carbonyl]amino]-benzoic acid
MF: C₂₁H₂₁NO₄
FW: 351.4
Purity: ≥98%
UV/Vis.: λ_{max}: 289 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

BMS 753 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 753 in the solvent of choice, which should be purged with an inert gas. BMS 753 is soluble in organic solvents such as ethanol and DMSO.

Description

BMS 753 is a synthetic retinoid and retinoic acid receptor α (RAR α) agonist ($K_i = 2$ nM).¹ It is selective for RAR α over RAR γ at 1 μ M. BMS 753 (10 and 100 nM) induces morphological differentiation of P19 embryonal carcinoma cells, as well as F9 embryonal carcinoma cells when used at a concentration of 100 nM in combination with the pan-retinoid X receptor (RXR) agonist BMS649.²

References

1. Géhin, M., Vivat, V., Wurtz, J.M., *et al.* Structural basis for engineering of retinoic acid receptor isotype-selective agonists and antagonists. *Chem. Biol.* **6(8)**, 519-529 (1999).
2. Taneja, R., Roy, B., Plassat, J.L., *et al.* Cell-type and promoter-context dependent retinoic acid receptor (RAR) redundancies for RAR β 2 and Hoxa-1 activation in F9 and P19 cells can be artefactually generated by gene knockouts. *Proc. Natl. Acad. Sci. USA* **93(12)**, 6197-6202 (1996).

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, 10/18/2022

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