

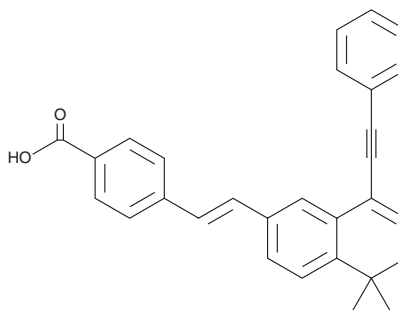
# PRODUCT INFORMATION



## BMS 493

Item No. 17418

**CAS Registry No.:** 215030-90-3  
**Formal Name:** 4-[(1E)-2-[5,6-dihydro-5,5-dimethyl-8-(2-phenylethynyl)-2-naphthalenyl]ethenyl]-benzoic acid  
**Synonym:** BMS 204493  
**MF:** C<sub>29</sub>H<sub>24</sub>O<sub>2</sub>  
**FW:** 404.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 239, 297, 328 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 493 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 493 in the solvent of choice. BMS 493 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of BMS 493 in these solvents is approximately 25 and 30 mg/ml, respectively.

BMS 493 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BMS 493 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BMS 493 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Nuclear retinoic acid receptors (RARs) are transcriptional regulators with roles in cell proliferation and differentiation.<sup>1,2</sup> BMS 493 is a pan-RAR inverse agonist that blocks RARα activity with an IC<sub>50</sub> value of 114 nM.<sup>3,4</sup> In all RAR types (RARα, RARβ, and RARγ), BMS 493 prevents the recruitment of transcriptional coactivators to RARs while stabilizing corepressor interactions.<sup>5,6</sup> BMS 493 has been used to elucidate the critical roles of RARs in development and immune response.<sup>3,7-9</sup>

### References

1. Duong, V. and Rochette-Egly, C. *Biochim. Biophys. Acta* **1812**(8), 1023-1031 (2011).
2. Rochette-Egly, C. and Germain, P. *Nucl. Recept. Signal.* **7**, 1-18 (2009).
3. Chazaud, C., Chambon, P., and Dollé, P. *Development* **126**(12), 2589-2596 (1999).
4. Garçça-Rodríguez, J., Pérez-Rodríguez, S., Ortiz, M.A., et al. *Bioor. Med. Chem.* **22**(4), 1285-1302 (2014).
5. le Maire, A., Teyssier, C., Erb, C., et al. *Nat. Struct. Mol. Biol.* **17**(7), 801-807 (2010).
6. Germain, P., Gaudon, C., Pogenberg, V., et al. *Chem. Biol.* **16**(5), 479-489 (2009).
7. Mollard, R., Ghyselinck, N.B., Wendling, O., et al. *Int. J. Dev. Biol.* **44**(5), 457-462 (2000).
8. Chazaud, C., Dollé, P., Rossant, J., et al. *Mech. Dev.* **120**(6), 691-700 (2003).
9. Geissmann, F., Revy, P., Brousse, N., et al. *J. Exp. Med.* **198**(4), 623-634 (2003).

#### 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, 02/14/2024

#### 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