

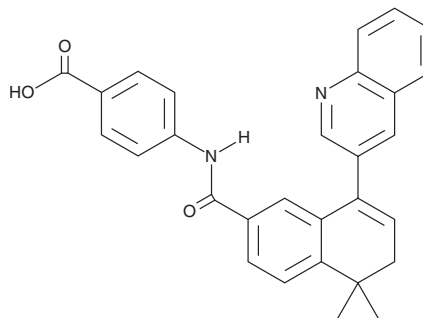
# PRODUCT INFORMATION



## BMS 195614

Item No. 16029

**CAS Registry No.:** 182135-66-6  
**Formal Name:** 4-[[[5,6-dihydro-5,5-dimethyl-8-(3-quinolinyl)-2-naphthalenyl]carbonyl]amino]-benzoic acid  
**Synonym:** BMS 614  
**MF:** C<sub>29</sub>H<sub>24</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 448.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 211, 236, 282 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 195614 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMS 195614 in the solvent of choice, which should be purged with an inert gas. BMS 195614 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BMS 195614 in these solvents is approximately 30 mg/ml.

BMS 195614 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BMS 195614 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BMS 195614 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

BMS 195614 is a neutral retinoic acid receptor (RAR)  $\alpha$ -selective antagonist ( $K_i = 2.5$  nM).<sup>1</sup> It antagonizes agonist-induced coactivator recruitment and moderately decreases SMRT binding to RAR but does not significantly affect nuclear receptor corepressor binding.<sup>2</sup>

### 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. Germain, P., Gaudon, C., Pogenberg, V., *et al.* Differential action on coregulator interaction defines inverse retinoid agonists and neutral antagonists. *Chem. Biol.* **16**(5), 479-489 (2009).

#### 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, 11/16/2022

#### 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