

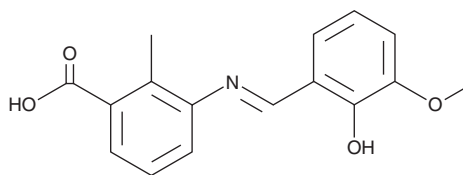
PRODUCT INFORMATION



C29

Item No. 27029

CAS Registry No.: 363600-92-4
Formal Name: 3-[[[(2-hydroxy-3-methoxyphenyl)methylene]amino]-2-methylbenzoic acid
Synonym: TLR2-IN-C29
MF: C₁₆H₁₅NO₄
FW: 285.3
Purity: ≥95%
UV/Vis.: λ_{max}: 226, 277, 320 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

C29 is supplied as a crystalline solid. A stock solution may be made by dissolving the C29 in the solvent of choice, which should be purged with an inert gas. C29 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of C29 in these solvents is approximately 0.5, 11, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of C29 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of C29 in PBS (pH 7.2) is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

C29 is an inhibitor of toll-like receptor 2 (TLR2) signaling.¹ It inhibits synthetic and bacterial TLR2 agonist-induced TLR2/1 and TLR2/6 signaling in HEK-293T cells expressing human TLR2 (HEK-TLR2) and THP-1 cells when used at a concentration of 50 μM. C29 also inhibits TLR2/1 signaling in primary murine macrophages without inducing cytotoxicity. It decreases the expression of IL-8 induced by the synthetic bacterial lipopeptide TLR2 agonist Pam₃CSK₄ (Item No. 24126) in HEK-TLR2 cells. C29 (150 μM) inhibits the Pam₃CSK₄-induced interaction between TLR2 and the adapter protein MyD88, blocks MAPK activation, and decreases NF-κB activation and IκBα degradation in THP-1 cells.

Reference

1. Mistry, P., Laird, M.H., Schwarz, R.S., *et al.* Inhibition of TLR2 signaling by small molecule inhibitors targeting a pocket within the TLR2 TIR domain. *Proc. Natl. Acad. Sci. U.S.A.* **112**(17), 5455-5460 (2015).

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, 12/19/2022

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