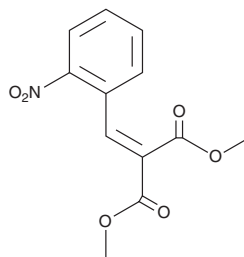


# PRODUCT INFORMATION



**NCI 126224**  
Item No. 39520

**CAS Registry No.:** 65974-52-9  
**Formal Name:** 2-[(2-nitrophenyl)methylene]-propanedioic acid, 1,3-dimethyl ester  
**Synonym:** NSC 126224  
**MF:** C<sub>12</sub>H<sub>11</sub>NO<sub>6</sub>  
**FW:** 265.2  
**Purity:** ≥98%  
**Supplied as:** A 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

NCI 126224 is supplied as a solid. A stock solution may be made by dissolving the NCI 126224 in the solvent of choice, which should be purged with an inert gas. NCI 126224 is soluble in chloroform and methanol.

## Description

NCI 126224 is a toll-like receptor 4 (TLR4) antagonist.<sup>1</sup> NCI 126224 selectively inhibits nitric oxide (NO) production induced by the TLR4 agonist LPS (IC<sub>50</sub> = 0.31 μM) over the TLR7/8 agonist R-848 (Item No. 14806), TLR1/2 agonist Pam<sub>3</sub>CSK<sub>4</sub> (Item No. 24126), and TLR3 agonist poly(I:C) in RAW 264.7 macrophages but also inhibits NO production induced by the TLR2/6 agonist FSL-1 (Item No. 24220) in the same cells at 0.6 μM. NCI 126224 inhibits LPS-induced NF-κB activity in a reporter assay using BV-2 microglial cells and LPS-induced increases in IL-1β and TNF-α levels in RAW 264.7 macrophages (IC<sub>50</sub>s = 5.92, 0.42, and 1.54 μM, respectively).

## Reference

1. Zhang, S., Cheng, K., Wang, X., *et al.* Selection, synthesis, and anti-inflammatory evaluation of the arylidene malonate derivatives as TLR4 signaling inhibitors. *Bioorg. Med. Chem.* **20(20)**, 6073-6079 (2012).

**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, 10/16/2023

## 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