

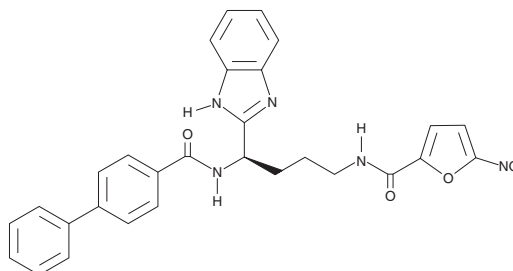
# PRODUCT INFORMATION



## LB244

Item No. 39832

**Formal Name:** (R)-N-(4-([1,1'-biphenyl]-4-carboxamido)-4-(1H-benzo[d]imidazol-2-yl)butyl)-5-nitrofuran-2-carboxamide  
**MF:** C<sub>29</sub>H<sub>25</sub>N<sub>5</sub>O<sub>5</sub>  
**FW:** 523.5  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥3 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

LB244 is supplied as a solid. A stock solution may be made by dissolving the LB244 in the solvent of choice, which should be purged with an inert gas. LB244 is soluble in chloroform.

### Description

LB244 is an irreversible STING antagonist.<sup>1</sup> It binds to STING in a cell-free assay and inhibits STING activation induced by the STING agonist diABZI in cells (EC<sub>50</sub> = 0.8 nM). LB244 (1 μM) inhibits diABZI-induced STING oligomerization and phosphorylation of IFN regulatory factor 3 (IRF3) and TANK-binding kinase 1 (TBK1) in primary mouse bone marrow-derived macrophages (BMDMs). *In vivo*, LB244 (5 mg/kg, i.p.) reduces diABZI-induced increases in serum IL-6 and IFN-β levels in mice.

### Reference

1. Barasa, L., Chaudhuri, S., Zhou, J.Y., *et al.* Development of LB244, an irreversible STING antagonist. *J. Am. Chem. Soc.* **145**(37), 20273-20288 (2023).

**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, 01/19/2024

### 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