

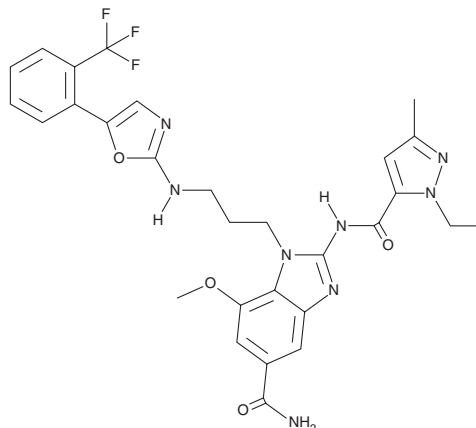
PRODUCT INFORMATION



STING Agonist D61

Item No. 38706

CAS Registry No.: 2850251-27-1
Formal Name: 2-[[[(1-ethyl-3-methyl-1H-pyrazol-5-yl)carbonyl]amino]-7-methoxy-1-[3-fluoromethyl]phenyl]-2-oxazolyl]amino]propyl]
Synonyms: D61, Stimulator of Interferon Genes Agonist D61
MF: C₂₉H₂₉F₃N₈O₄
FW: 610.6
Purity: ≥95%
UV/Vis.: λ_{max}: 322 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

STING agonist D61 is supplied as a solid. A stock solution may be made by dissolving the STING agonist D61 in the solvent of choice, which should be purged with an inert gas. STING agonist D61 is soluble in methanol and DMSO.

Description

STING agonist D61 (D61) is an agonist of stimulator of interferon genes (STING).¹ It induces expression of an IFN3-inducible secreted alkaline phosphatase (SEAP) reporter gene and an IFN-β-inducible reporter gene in cell-based assays (EC₅₀s = 52.9 and 116 nM, respectively). D61 (4, 6, and 8 μM) increases the expression of mRNA encoding IFN-β and chemokine (C-X-C motif) ligand 10 (CXCL10) and phosphorylation of TANK-binding kinase 1 (TBK1), IRF3, and STING in THP-1 monocytes. *In vivo*, D61 (0.25 mg/kg every other day) reduces tumor volume without affecting body weight in a CT26 murine colon cancer model.

Reference

1. Chen, N.-N., Zhang, H., Zhu, Q.-S., *et al.* Development of orally bioavailable amidobenzimidazole analogues targeting stimulator of interferon gene (STING) receptor. *J. Med. Chem.* **66**(8), 5584-5610 (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, 10/17/2023

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