

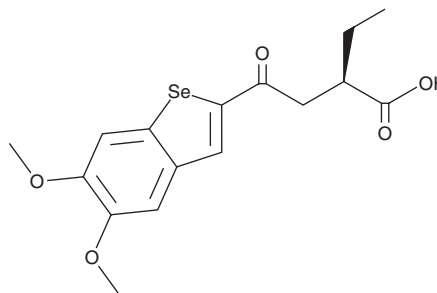
PRODUCT INFORMATION



BSP-16

Item No. 37992

CAS Registry No.: 2727249-47-8
Formal Name: αR-ethyl-5,6-dimethoxy-γ-oxo-benzo[b]selenophene-2-butanoic acid
Synonym: LF250
MF: C₁₆H₁₈O₅Se
FW: 369.3
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 332 nm
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.

Description

BSP-16 is a stimulator of interferon genes (STING) agonist.¹ It binds to STING ($K_d = 23 \mu\text{M}$) and induces the production of IFN- β , IL-6, and chemokine (C-X-C motif) ligand 10 (CXCL10) in THP-1 cells when used at concentrations of 10, 25, and 50 μM . BSP-16 (50 μM) induces phosphorylation of IFN regulatory factor 3 (IRF3) and TANK-binding kinase 1 (TBK1) in THP-1 cells. *In vivo*, BSP-16 (15 and 30 mg/kg) reduces tumor volume in an MC-38 murine carcinoma model and induces tumor regression in a CT26 murine colorectal carcinoma model. It prevents tumor cell re-challenge-induced tumor formation in a CT26 murine carcinoma model. BSP-16 (20 mg/kg) also reduces tumor volume and increases the number of activated CD8⁺ T cells and natural killer (NK) cells in a 4T1 murine mammary carcinoma model.

Reference

1. Feng, X., Pan, L., Qian, Z., *et al.* Discovery of selenium-containing STING agonists as orally available antitumor agents. *J. Med. Chem.* **65**(22), 15048-15065 (2022).

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, 06/16/2023

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