



Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

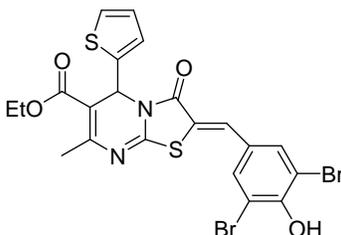
Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Product Name: **IspF Inhibitor 1**
(2Z)-2-(3,5-Dibromo-4-hydroxybenzylidene)-7-methyl-3-oxo-5-(2-thienyl)-2,3-dihydro-5H-[1,3]thiazolo[3,2-a]pyrimidine-6-carboxylate

Product Number: **B-4102**



Molecular Formula: C₂₁H₁₆Br₂N₂O₄S₂ **MW:** 584.30 **CAS:** 1240789-72-3

Solubility: DMSO, CHCl₃:MeOH (9:1), >5 mg/mL

Storage: 4 °C

Field of Interest: The non-mevalonate, or MEP, pathway for isoprenoid biosynthesis is essential in gram-negative, some gram-positive bacteria, plants, and protozoa making it an attractive pathway to target for antibiotics since mammals use the mevalonate pathway to synthesize isoprenoids. IspF is the fourth step in the pathway and converts diphosphocytidyl-2-methylerythritol 2-phosphate into 2-methylerythritol 2,4-cyclodiphosphate (cMEPP). IspF inhibitor 1 (compound 3 in Geise *et al.*) has low micromolar activity (IC₅₀ = 6.1-32 μM) against the recombinant enzymes and inhibits the growth of *P. falciparum* in a red blood cell assay.

References: J.G. Geist *et al.* "Thiazolopyrimidine Inhibitors of 2-Methylerythritol 2,4-Cyclodiphosphate Synthase (IspF) from *Mycobacterium tuberculosis* and *Plasmodium falciparum*" *ChemMedChem* 2010, 5, 1092-1101.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. IspF Inhibitor 1 is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.