

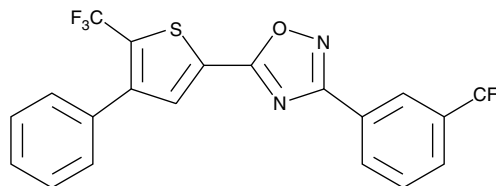
Product Information



SEW2871

Item No. 10006440

CAS Registry No.: 256414-75-2
Formal Name: 5-[4-phenyl-5-(trifluoromethyl)-2-thienyl]-3-[3-(trifluoromethyl)phenyl]-1,2,4-oxadiazole
MF: C₂₀H₁₀F₆N₂OS
FW: 440.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 242, 273 nm



Laboratory Procedures

For long term storage, we suggest that SEW2871 be stored as supplied at -20°C. It should be stable for at least two years.

SEW2871 is supplied as a crystalline solid. A stock solution may be made by dissolving the SEW2871 in an organic solvent purged with an inert gas. SEW2871 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SEW2871 in these solvents is 5, 2, and 30 mg/ml, respectively.

SEW2871 is sparingly soluble in aqueous buffers. Therefore, further dilutions of the organic solvent solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Sphingosine-1-phosphate receptor 1 (S1P₁) is one of five high affinity G protein-coupled S1P receptors which mediate a variety of effects including lymphocyte recirculation in the blood.¹ Non-selective S1P receptor agonists, such as FTY720, produce clinical immunosuppression useful for preventing transplant rejection and treating autoimmune diseases. However, they also cause bradycardia by activating S1P₃, the receptor responsible for regulation of heart rate.² SEW2871 is a selective S1P₁ receptor agonist in both human and mouse that is not active at the S1P₂₋₅ receptors. SEW2871, therefore, suppresses the immune response by decreasing the number of lymphocytes circulating in blood without causing bradycardia.²

References

1. Kluk, M.J. and Hla, T. Signaling of sphingosine-1-phosphate *via* the S1P/EDG-family of G-protein-coupled receptors. *Biochim. Biophys. Acta* **1582**, 72-80 (2002).
2. Sanna, M.G., Liao, J., Jo, E., *et al.* Sphingosine 1-phosphate (S1P) receptor subtypes SIP₁ and SIP₃, respectively, regulate lymphocyte recirculation and heart rate. *J. Biol. Chem.* **279**(14), 13839-13848 (2004).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10006440

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 06/18/2012

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com