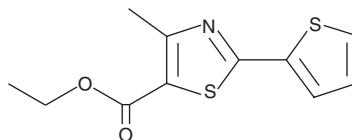


PRODUCT INFORMATION

ethyl 4-methyl-2-(2-thienyl)-thiazole-5-Carboxylate

Item No. 11056

CAS Registry No.: 56421-62-6
Formal Name: 4-methyl-2-(2-thienyl)-5-thiazolecarboxylic acid, ethyl ester
MF: C₁₁H₁₁NO₂S₂
FW: 253.3
Purity: ≥97%
UV/Vis.: λ_{max}: 239, 335 nm
Supplied as: A crystalline 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

ethyl 4-methyl-2-(2-thienyl)-thiazole-5-Carboxylate is supplied as a crystalline solid. A stock solution may be made by dissolving the ethyl 4-methyl-2-(2-thienyl)-thiazole-5-carboxylate in an organic solvent purged with an inert gas. ethyl 4-methyl-2-(2-thienyl)-thiazole-5-Carboxylate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ethyl 4-methyl-2-(2-thienyl)-thiazole-5-carboxylate in these solvents is approximately 1, 5, and 10 mg/ml, respectively.

ethyl 4-methyl-2-(2-thienyl)-thiazole-5-Carboxylate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ethyl 4-methyl-2-(2-thienyl)-thiazole-5-carboxylate should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ethyl 4-methyl-2-(2-thienyl)-thiazole-5-Carboxylate has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ethyl 4-methyl-2-(2-thienyl)-thiazole-5-Carboxylate is a synthetic intermediate useful for pharmaceutical synthesis.

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, 11/09/2017

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