

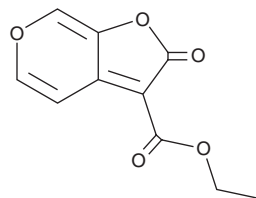
PRODUCT INFORMATION



Karrikinolide 3-ethyl ester

Item No. 31675

CAS Registry No.: 1030604-40-0
Formal Name: 2-oxo-2H-furo[2,3-c]pyran-3-carboxylic acid, ethyl ester
MF: $C_{10}H_8O_5$
FW: 208.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

Karrikinolide 3-ethyl ester is supplied as a solid. A stock solution may be made by dissolving the karrikinolide 3-ethyl ester in the solvent of choice, which should be purged with an inert gas. Karrikinolide 3-ethyl ester is soluble in methanol and DMSO.

Description

Karrikinolide 3-ethyl ester is a synthetic intermediate.¹ It has been used in the synthesis of butanolide seed germination stimulants.

Reference

1. Sun, K., Chen, Y., Wagerle, T., *et al.* Synthesis of butenolides as seed germination stimulants. *Tetrahedron Lett.* **49**, 2922-2925 (2008).

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, 11/10/2022

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