

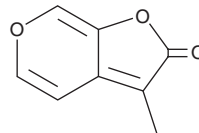
# PRODUCT INFORMATION



## Karrikinolide

Item No. 29935

CAS Registry No.: 857054-02-5  
Formal Name: 3-methyl-2H-furo[2,3-c]pyran-2-one  
Synonym: Karrikin 1  
MF: C<sub>8</sub>H<sub>6</sub>O<sub>3</sub>  
FW: 150.1  
Purity: ≥98%  
Supplied as: A 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

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

### Description

Karrikinolide is a plant growth regulator that has been found in plant-derived smoke.<sup>1</sup> It increases the germination rate and seedling mass in a panel of eight arable weed species when applied to filter paper at a concentration of 1 μM. Topical application of karrikinolide (0.1 nM) increases the number of leaves, leaf length, and fresh and dry leaf weights in commercial onion (*A. cepa*) plants.<sup>2</sup>

### References

1. Daws, M.I., Davies, J., Pritchard, H.W., *et al.* Butenolide from plant-derived smoke enhances germination and seedling growth of arable weed species. *Plant Growth Regul.* **51**, 73-82 (2006).
2. Kulkarni, M.G., Ascough, G.D., Verschaeve, L., *et al.* Effect of smoke-water and a smoke-isolated butenolide on the growth and genotoxicity of commercial onion. *Sci. Hortic. (Amsterdam)* **124(4)**, 434-439 (2010).

#### 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, 03/25/2020

#### 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