

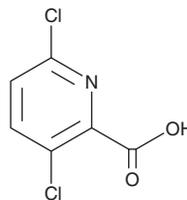
PRODUCT INFORMATION



Clopyralid

Item No. 24229

CAS Registry No.: 1702-17-6
Formal Name: 3,6-dichloro-2-pyridinecarboxylic acid
MF: C₆H₃Cl₂NO₂
FW: 192.0
Purity: ≥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

Clopyralid is supplied as a solid. A stock solution may be made by dissolving the clopyralid in the solvent of choice, which should be purged with an inert gas. Clopyralid is slightly soluble in DMSO.

Description

Clopyralid is an herbicide that selectively inhibits the growth of the broadleaf weed *L. palustris* over a panel of 17 New Zealand native non-broadleaf plant species when applied at a concentration of 1.5 mg/L.¹ It does not inhibit growth of human pluripotent stem cells (IC₅₀ = >100 μM) or decrease nuclear translocation of the transcription factor SOX17, a marker of teratogenic risk.² Formulations containing clopyralid have been used as herbicides in agriculture.

References

1. Champion, P.D. Selective control of weeds in New Zealand wetlands. *Proceedings of the 51st New Zealand Plant Protection Conference* 251-254 (1998).
2. Kameoka, S., Babiarz, J., Kolaja, K., *et al.* A high-throughput screen for teratogens using human pluripotent stem cells. *Toxicol. Sci.* **137**(1), 76-90 (2014).

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, 09/19/2023

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