

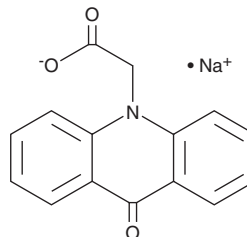
PRODUCT INFORMATION



Cridanimod (sodium salt)

Item No. 27884

CAS Registry No.: 58880-43-6
Formal Name: 9-oxo-10(9H)-acridineacetic acid, monosodium salt
Synonym: 10-Carboxymethyl-9-Acridanone
MF: C₁₅H₁₀NO₃ • Na
FW: 275.2
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 255, 400 nm
Supplied as: A crystalline 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

Cridanimod (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the cridanimod (sodium salt) in the solvent of choice, which should be purged with an inert gas. Cridanimod (sodium salt) is soluble in the organic solvent DMSO, at a concentration of approximately 20 mg/ml.

Description

Cridanimod is an inducer of type I interferon (IFN) production.^{1,2} It induces IRF3 phosphorylation, IFN-β production, and NF-κB activation in wild-type, but not in stimulator of interferon genes (STING) mutant, murine macrophages.¹ *In vivo*, cridanimod (112-1,792 mg/kg) increases plasma levels of IFN in weanling and adult mice.² Cridanimod inhibits viral infection in mouse models of Semliki forest, coxsackie B1, Columbia SK, herpes, and pseudorabies viruses with protective doses (PD₅₀s) ranging from 17-320 mg/kg.³ It increases uterine expression of estrogen and progesterone receptors in ovariectomized rats.⁴ Cridanimod also reverses tamoxifen-induced decreases in progesterone receptor expression in young rats.

References

1. Cavlar, T., Deimling, T., Ablasser, A., *et al.* Species-specific detection of the antiviral small-molecule compound CMA by STING. *EMBO J.* **32(10)**, 1440-1450 (2013).
2. Taylor, J.L., Schoenherr, C.K., and Grossberg, S.E. High-yield interferon induction by 10-carboxymethyl-9-acridanone in mice and hamsters. *Antimicrob. Agents Chemother.* **18(1)**, 20-26 (1980).
3. Kramer, M.J., Cleeland, R., and Grunberg, E. Antiviral activity of 10-carboxymethyl-9-acridanone. *Antimicrob. Agents Chemother.* **9(2)**, 233-238 (1976).
4. Surkov, K.G., Tsyrlina, E.V., Konstantinova, M.M., *et al.* Neovir, an interferon inducer, modifies expression of steroid hormone receptors in hormone-dependent tissues and restores sensitivity to tamoxifen in patients with inoperable breast cancer. *Vopr. Onkol.* **42(6)**, 28-32 (1996).

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, 12/07/2022

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