

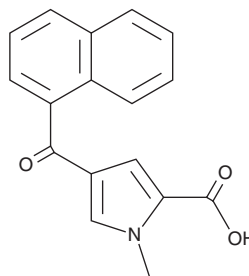
PRODUCT INFORMATION



Napirimus

Item No. 41873

CAS Registry No.: 70696-66-1
Formal Name: 1-methyl-4-(1-naphthalenylcarbonyl)-1H-pyrrole-2-carboxylic acid
MF: $C_{17}H_{13}NO_3$
FW: 279.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Napirimus is supplied as a solid. A stock solution may be made by dissolving the napirimus in the solvent of choice, which should be purged with an inert gas. Napirimus is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Napirimus is an immunosuppressant.¹ It is efficacious and preventative in a rat model of polyarthritis induced by complete Freund's adjuvant when administered at doses of 200 and 100 mg/kg, respectively. Napirimus is 50 and 90% protective against experimental autoimmune encephalomyelitis (EAE), previously known as allergic encephalomyelitis, in rats when administered at doses of 100 and 400 mg/kg, respectively.

Reference

1. Bessin, P., Mazarin, C., and Bonnet, J. Pharmaceutical compositions having immuno-suppressive properties. *Albert Rolland, S.A. 4,617,315*, (1986).

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, 08/29/2024

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