

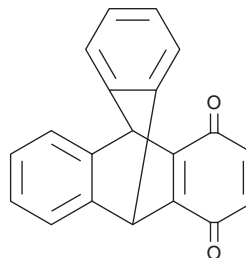
PRODUCT INFORMATION



NFAT Activation Inhibitor III

Item No. 21812

CAS Registry No.: 3519-82-2
Formal Name: 9,10-dihydro-9,10[1',2']-benzenoanthracene-1,4-dione
Synonyms: INCA-6, NSC 25996, Nuclear Factor of Activated T cells Inhibitor III
MF: C₂₀H₁₂O₂
FW: 284.3
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 252 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

NFAT activation inhibitor III is supplied as a crystalline solid. A stock solution may be made by dissolving the NFAT activation inhibitor III in the solvent of choice, which should be purged with an inert gas. NFAT activation inhibitor III is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of NFAT activation inhibitor III in these solvents is approximately 2 and 20 mg/ml, respectively.

Description

NFAT activation inhibitor III is a cell-permeant inhibitor of signaling through nuclear factor of activated T cells (NFAT).¹ It binds the substrate recognition site of the phosphatase calcineurin ($K_d = 0.8 \mu\text{M}$), preventing it from binding and dephosphorylating NFAT.¹ NFAT activation inhibitor III blocks the nuclear import of NFAT in T cells stimulated with ionomycin (Item No. 10004974). It prevents the induced expression of cytokine mRNAs in T cells stimulated with phorbol ester and ionomycin.¹

Reference

1. Roehrl, M.H.A., Kang, S., Aramburu, J., *et al.* Selective inhibition of calcineurin-NFAT signaling by blocking protein-protein interaction with small organic molecules. *Proc. Natl. Acad. Sci. USA* **101**(20), 7554-7559 (2004).

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, 12/08/2022

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