

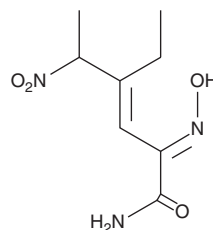
PRODUCT INFORMATION



FK-409

Item No. 82290

CAS Registry No.: 138472-01-2
Formal Name: 4-ethyl-2E-(hydroxyimino)-5-nitro-3E-hexenamamide
Synonyms: NOR-3
MF: C₈H₁₃N₃O₄
FW: 215.2
Purity: ≥98%
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

FK-409 is supplied as a crystalline solid. A stock solution may be made by dissolving the FK-409 in the solvent of choice, which should be purged with an inert gas. FK-409 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of FK-409 in these solvents is approximately 36, 20, and 34 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of FK-409 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of FK-409 in PBS (pH 7.2) is approximately 2.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

FK-409 is a NO donor. It liberates NO in a time-dependent manner with a half-life of approximately 40 minutes in phosphate buffer (pH 7.4) at 37°C.¹ The IC₅₀ for the inhibition of platelet aggregation is 4.9 μM.²

References

1. Kita, Y., Hirasawa, Y., Maeda, K., *et al.* Spontaneous nitric oxide release accounts for the potent pharmacological actions of FK409. *Eur. J. Pharmacol.* **257(1-2)**, 123-130 (1994).
2. Fukuyama, S., Hirasawa, Y., Kato, Y., *et al.* Structure-activity relationships of spontaneous nitric oxide releasers, FK409 and its derivatives. *J. Pharmacol. Exp. Ther.* **282(1)**, 236-242 (1997).

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, 11/03/2022

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