

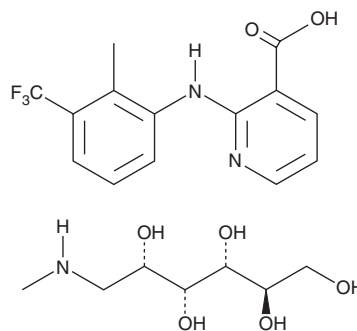
PRODUCT INFORMATION



Flunixin (meglumine)

Item No. 26644

CAS Registry No.: 42461-84-7
Formal Name: 1-deoxy-1-(methylamino)-D-glucitol, 2-[[2-methyl-3-(trifluoromethyl)phenyl]amino]-3-pyridinecarboxylate
Synonym: NIH 10250
MF: C₁₄H₁₁F₃N₂O₂ • C₇H₁₇NO₅
FW: 491.5
Purity: ≥98%
UV/Vis.: λ_{max}: 279 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

Flunixin (meglumine) is supplied as a crystalline solid. A stock solution may be made by dissolving the flunixin (meglumine) in the solvent of choice. Flunixin (meglumine) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of flunixin (meglumine) in ethanol is approximately 0.25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

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 flunixin (meglumine) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of flunixin (meglumine) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Flunixin is a non-steroidal anti-inflammatory drug (NSAID) that inhibits COX-1 and COX-2 (IC₅₀s = 3.24 and 0.55 μM, respectively).¹ It inhibits carrageenan-induced prostaglandin E₂ (PGE₂; Item No. 14010) and thromboxane B₂ (TXB₂; Item No. 19030) production in sheep exudate and serum, respectively, when administered at a dose of 1.1 mg/kg.² Flunixin (1.1 mg/l/kg) has anti-inflammatory, analgesic, and antipyretic activity in horses.³ Formulations containing flunixin have been used in the veterinary treatment of pain and fever in livestock.

References

1. Bryant, C.E., Farnfield, B.A., and Janicke, H.J. Evaluation of the ability of carprofen and flunixin meglumine to inhibit activation of nuclear factor kappa B. *Am. J. Vet. Res.* **64**(2), 211-215 (2003).
2. Cheng, Z., Nolan, A.M., and McKellar, Q.A. Measurement of cyclooxygenase inhibition in vivo: A study of two non-steroidal anti-inflammatory drugs in sheep. *Inflammation* **22**(4), 353-366 (1998).
3. Houdeshell, J.W. and Hennessey, P.W. A new nonsteroidal, anti-inflammatory analgesic for horses. *J. Equine Vet Sci.* **1**(2), 57-63 (1977).

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, 12/19/2022

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