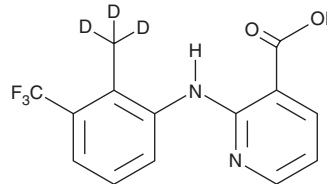


PRODUCT INFORMATION



Flunixin-d₃ Item No. 35056

CAS Registry No.: 1015856-60-6
Formal Name: 2-[[2-(methyl-d₃)-3-(trifluoromethyl)phenyl]amino]-3-pyridinecarboxylic acid
MF: C₁₄H₈D₃F₃N₂O₂
FW: 299.3
Chemical Purity: ≥90% (Flunixin)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A 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-d₃ is intended for use as an internal standard for the quantification of flunixin (Item No. 26644) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Flunixin-d₃ is supplied as a solid. A stock solution may be made by dissolving the flunixin-d₃ in the solvent of choice, which should be purged with an inert gas. Flunixin-d₃ is soluble in DMSO.

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 per 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, 10/17/2022

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