

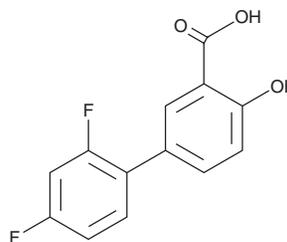
PRODUCT INFORMATION



Diflunisal

Item No. 19187

CAS Registry No.: 22494-42-4
Formal Name: 2',4'-difluoro-4-hydroxy-[1,1'-biphenyl]-3-carboxylic acid
Synonyms: 5-(2,4-Difluorophenyl)salicylic Acid, MK-647
MF: C₁₃H₈F₂O₃
FW: 250.2
Purity: ≥98%
UV/Vis.: λ_{max}: 228, 253 nm
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥4 years



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

Laboratory Procedures

Diflunisal is supplied as a crystalline solid. A stock solution may be made by dissolving the diflunisal in the solvent of choice. Diflunisal is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of diflunisal in ethanol and DMSO is approximately 25 mg/ml and approximately 30 mg/ml in DMF.

Diflunisal is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, diflunisal should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Diflunisal has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Diflunisal is a non-steroidal anti-inflammatory drug (NSAID) that inhibits both COX-1 (IC₅₀ = 113 μM) and COX-2 (IC₅₀s = 8.2 and 134 μM for human whole blood assay and human-modified whole blood assays, respectively).¹ Peak plasma levels are achieved within two hours, with little metabolism before excretion in the urine.² The terminal plasma half-life is approximately eight hours.²

References

1. Warner, T.D., Giuliano, F., Vojnovic, I., *et al.* Nonsteroid drug selectivities for cyclo-oxygenase-1 rather than cyclo-oxygenase-2 are associated with human gastrointestinal toxicity: A full *in vitro* analysis. *Proc. Natl. Acad. Sci. USA* **96**, 7563-7568 (1999).
2. Tempero, K.F., Cirillo, V.J., and Steelman, S.L. Diflunisal: A review of pharmacokinetic and pharmacodynamic properties, drug interactions, and special tolerability studies in humans. *Br. J. Clin. Pharmacol.* **4 Suppl 1**, 31S-36S (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, 11/11/2022

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