

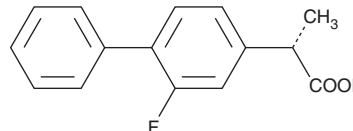
PRODUCT INFORMATION



(S)-Flurbiprofen

Item No. 10004207

CAS Registry No.: 51543-39-6
Formal Name: (S)-(+)-2-fluoro- α -methyl-4-biphenylacetic acid
MF: C₁₅H₁₃FO₂
FW: 244.3
Purity: $\geq 99\%$
UV/Vis.: λ_{max} : 247 nm
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥ 2 years



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

Laboratory Procedures

(S)-Flurbiprofen is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-flurbiprofen in an organic solvent purged with an inert gas. (S)-Flurbiprofen is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (S)-flurbiprofen in these solvents is at least 10 mg/ml in DMSO and 25 mg/ml in ethanol 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 (S)-flurbiprofen can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of (S)-flurbiprofen in PBS (pH 7.2) is at least 0.9 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(S)-Flurbiprofen is the COX-active enantiomer of the non-selective COX inhibitor flurbiprofen (Item No. 70250) with IC₅₀ values of 0.48 and 0.47 μM for COX-1 and COX-2, respectively, in guinea pig whole blood.¹ It inhibits the release of 6-keto prostaglandin F_{1 α} (6-keto-PGF_{1 α} ; Item No. 15210) and thromboxane B₂ (TXB₂; Item No. 19030) from rat whole blood, gastric mucosa, lung, and jejunal tissue *ex vivo* in a dose-dependent manner.² (S)-Flurbiprofen (1 nM) inhibits basal and bradykinin-, serotonin-, and histamine-stimulated prostaglandin E₂ (PGE₂) release from isolated skin flaps of rat lower hind paws.³ It also inhibits release of the neuroinflammatory marker calcitonin gene-related peptide (CGRP; Item Nos. 24405 | 24725 | 24728) when used at a concentration of 1 μM . *In vivo*, (S)-flurbiprofen reduces the number of flinches per minute in the formalin test in rats, indicating antinociceptive activity.⁴

References

1. Carabaza, A., Cabre, F., Rotllan, E., *et al.* Stereoselective inhibition of inducible cyclooxygenase by chiral nonsteroidal antiinflammatory drugs. *J. Clin. Pharmacol* **36**(6), 505-512 (1996).
2. Peskar, B.M., Kluge, S., Peskar, B.A., *et al.* Effects of pure enantiomers of flurbiprofen in comparison to racemic flurbiprofen on eicosanoid release from various rat organs *ex vivo*. *Prostaglandins* **42**(6), 515-531 (1991).
3. Averbeck, B., Peisler, M., Izydorczyk, I., *et al.* Inflammatory mediators do not stimulate CGRP release if prostaglandin synthesis is blocked by S(+)-flurbiprofen in isolated rat skin. *Inflamm. Res.* **52**(12), 519-523 (2003).
4. Malmberg, A.B. and Yaksh, T.L. Antinociception produced by spinal delivery of the S and R enantiomers of flurbiprofen in the formalin test. *Eur. J. Pharmacol.* **256**(2), 205-209 (1994).

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, 06/04/2018

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