

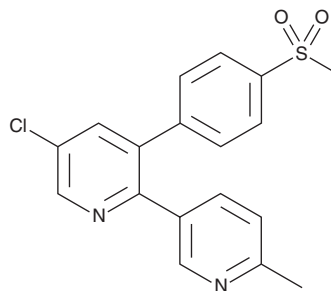
# PRODUCT INFORMATION



## Etoricoxib

Item No. 10091

**CAS Registry No.:** 202409-33-4  
**Formal Name:** 5-chloro-6'-methyl-3-[4-(methylsulfonyl)phenyl]-2,3'-bipyridine  
**Synonyms:** L-791,456, MK-0663  
**MF:** C<sub>18</sub>H<sub>15</sub>ClN<sub>2</sub>O<sub>2</sub>S  
**FW:** 358.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 236, 285 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Special Conditions:** Fatal in contact with skin



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

### Laboratory Procedures

Etoricoxib is supplied as a crystalline solid. A stock solution may be made by dissolving the etoricoxib in the solvent of choice. Etoricoxib is freely soluble in organic solvents such as methanol and DMSO, which should be purged with an inert gas.

### Description

Etoricoxib is a dipyrindinyl compound that demonstrates high *in vitro* and *ex vivo* selectivity for COX-2 over COX-1 in several assays, e.g., in the production of PGE<sub>2</sub> by CHO cells expressing either COX-2 (IC<sub>50</sub> = 79 nM) or COX-1 (IC<sub>50</sub> > 50 μM).<sup>1,2</sup> Oral etoricoxib is well absorbed and metabolized extensively via oxidation, with metabolites excreted largely in the urine.<sup>3</sup>

### References

1. Riendeau, D., Percival, M.D., Brideau, C., *et al.* Etoricoxib (MK-0663): Preclinical profile and comparison with other agents that selectively inhibit cyclooxygenase-2. *J. Pharmacol. Exp. Ther.* **296**, 558-566 (2001).
2. Dallob, A., Hawkey, C.J., Greenberg, H., *et al.* Characterization of etoricoxib, a novel, selective COX-2 inhibitor. *J. Clin. Pharmacol.* **43(6)**, 573-585 (2003).
3. Rodrigues, A.D., Halpin, R.A., Geer, L.A., *et al.* Absorption, metabolism, and excretion of etoricoxib, a potent and selective cyclooxygenase-2 inhibitor, in healthy male volunteers. *Drug Metab. Dispos.* **31(2)**, 224-232 (2003).

#### 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, 11/08/2022

#### 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