

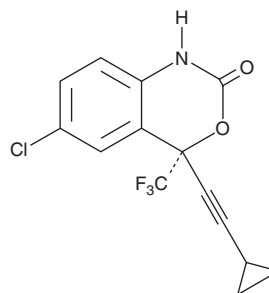
# PRODUCT INFORMATION



## Efavirenz

Item No. 14412

**CAS Registry No.:** 154598-52-4  
**Formal Name:** (4S)-6-chloro-4-(2-cyclopropylethynyl)-1,4-dihydro-4-(trifluoromethyl)-2H-3,1-benzoxazin-2-one  
**Synonyms:** DMP 266, EFV, L-743,726  
**MF:** C<sub>14</sub>H<sub>9</sub>ClF<sub>3</sub>NO<sub>2</sub>  
**FW:** 315.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 247, 294 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

Efavirenz is supplied as a crystalline solid. A stock solution may be made by dissolving the efavirenz in the solvent of choice, which should be purged with an inert gas. Efavirenz is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of efavirenz in ethanol and DMF is approximately 20 mg/ml and approximately 14 mg/ml in DMSO.

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

### Description

Efavirenz is a non-nucleoside reverse transcriptase inhibitor (NNRTI) that binds to wild-type and mutant HIV-1 RTs ( $K_i$ s = 2.93 and 3.85-56.5 nM, respectively).<sup>1</sup> It inhibits wild-type and mutant HIV-1 viral replication in MT-4 human T lymphoid cells ( $IC_{95}$ s = 1.5-1,500 nM). Efavirenz also prevents RNA plus-strand initiation with an  $IC_{50}$  value of 17 nM.<sup>2</sup> *In vivo*, efavirenz reduces HIV-1 cDNA in spleen of HIV-1-challenged HIV-susceptible transgenic rats.<sup>3</sup> Formulations containing efavirenz have been used in combination therapy for the treatment of HIV-1.<sup>4,5</sup>

### References

1. Young, S.D., Britcher, S.F., Tran, L.O., *et al.* L-743,726 (DMP-266): A novel, highly potent nonnucleoside inhibitor of the human immunodeficiency virus type 1 reverse transcriptase. *Antimicrob. Agents Chemother.* **39(12)**, 2602-2605 (1995).
2. Grobler, J.A., Dornadula, G., Rice, M.R., *et al.* HIV-1 reverse transcriptase plus-strand initiation exhibits preferential sensitivity to non-nucleoside reverse transcriptase inhibitors *in vitro*. *J. Biol. Chem.* **282(11)**, 8005-8010 (2007).
3. Goffinet, C., Allespach, I., and Keppler, O.T. HIV-susceptible transgenic rats allow rapid preclinical testing of antiviral compounds targeting virus entry or reverse transcription. *Proc. Natl. Acad. Sci. U.S.A.* **104(3)**, 1015-1020 (2007).
4. Sheran, M. The nonnucleoside reverse transcriptase inhibitors efavirenz and nevirapine in the treatment of HIV. *HIV Clin. Trials* **6(3)**, 158-168 (2005).
5. Rakhmanina, N.Y. and van den Anker, J.N. Efavirenz in the therapy of HIV infection. *Expert Opin. Drug Metab. Toxicol.* **6(1)**, 95-103 (2010).

#### 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