

# TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

---

## Etoposide

**Catalog #:** 4684-096-06

**Size:** 100 µl

**Description:** Etoposide complexes with topoisomerase II and DNA to enhance double-strand and single-strand cleavage of DNA and reversibly inhibit religation. Etoposide also blocks the cell cycle in S-phase and G2-phase; induces apoptosis in normal and tumor cell lines; inhibits synthesis of the oncoprotein Mdm2 and induces apoptosis in tumor lines that overexpress Mdm2.

**Formula weight:** 588.56

**Physical State:** Etoposide is provided in DMSO. The final reagent concentration is 10 mM.

**Storage:** Store at -20 °C.  
Bring to room temperature before use.

**Applications:** Induction of apoptosis *in vitro*. To induce apoptosis, dilute etoposide in media to a final concentration ranging from 10 µM to 100 µM. The concentration and length of treatment are cell type dependant and require optimization. Please refer to MSDS before handling this product.

## References:

1. Chow, K.C., and Ross, W.E., 1987. Topoisomerase-specific drug sensitivity in relation to cell cycle progression. *Mol. Cell Biol.*, **7**:3119-3123.
2. Fearnhead HO, M. Chwalinski , R.T. Snowden, M.G. Ormerod, GM.Cohen.1994. Dexamethasone and etoposide induce apoptosis in rat thymocytes from different phases of the cell cycle. *Biochem Pharmacol* **48**:1073-9.
3. Burden, D.A., et al., 1996. Topoisomerase II-etoposide interactions direct the formation of drug-induced enzyme-DNA cleavage complexes. *J. Biol. Chem.*, **271**:29238-29244.
4. Hande, K.R. 1998. Etoposide: four decades of development of a topoisomerase II inhibitor. *Eur.J. Cancer*, **34**:1514-1521.
5. Kaufmann, S.H., 1998. Cell death induced by topoisomerase-targeted drugs: more questions than answers. *Biochim. Biophys. Acta*, **1400**:195-211.

© 2018 Trevigen, Inc. All Rights Reserved. Trevigen is a registered trademark of Trevigen, Inc.

---

## TREVIGEN®

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: [info@trevigen.com](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

## **Etoposide**

**Catalog #: 4684-096-06**

**Storage: -20 °C**

**TREVIGEN®**

**1-800-873-8443**