

## SildENafil-D8 citRATE



### ORDERING INFORMATION

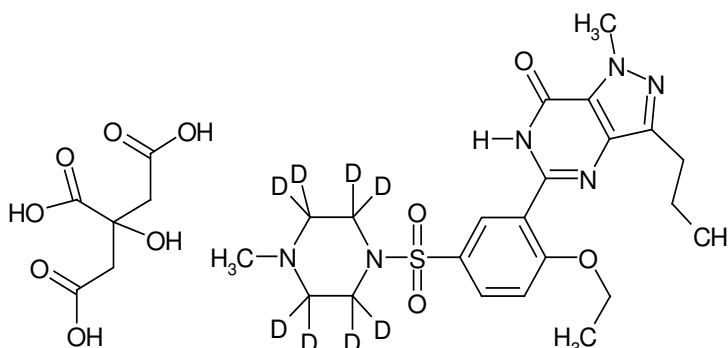
**Catalogue Number:** MI7948

**Size:** 1 mg  
5 mg  
10 mg

**Formulation:** solid

**Stability:** ≥ 1 year at +4 °C

**Storage:** +4 °C



$C_{28}H_{30}N_6O_{11}SD_8$

- **CAS NUMBER:** 171599-83-0 (for non labelled molecule)
- **MOLECULAR WEIGHT:** 674.8
- **PURITY:** ≥ 97% (%D: ≥ 98%)
- **APPLICATION:** Isotopically Labeled Pharmaceutical Reference Standard.
- **DESCRIPTION:**  
Sildenafil citrate is a drug used to treat male erectile dysfunction (impotence) and pulmonary arterial hypertension (PAH).
- **BIBLIOGRAPHY:**
  1. Orally active, selective type 5 cGMP phosphodiesterase inhibitor. Prepn: A. S. Bell *et al.*, EP 463756; *idem*, US 5250534 (1992, 1993 both to Pfizer).
  2. Structure-activity study: N. K. Terrett *et al.*, *Bioorg. Med. Chem. Lett.* 6, 1819 (1996). Clinical trial in impotence: I. Goldstein *et al.*, *N. Engl. J. Med.* 338, 1397 (1998).
  3. Review of pharmacology: M. Boolell *et al.*, *Int. J. Impot. Res.* 8, 47-52 (1996); of clinical experience: D. G. Hatzichristou, *ibid.* 14, Suppl. 1, S43-S52 (2002).
  4. Clinical evaluation in pulmonary hypertension: B. K. S. Sastry *et al.*, *J. Am. Coll. Cardiol.* 43, 1149 (2004).



February 2008

**Should you wish further information, do not hesitate to contact us.**