

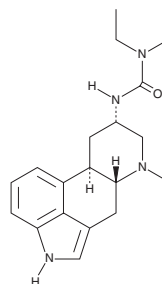
# PRODUCT INFORMATION



## Terguride

Item No. 42028

**CAS Registry No.:** 37686-84-3  
**Formal Name:** N,N-diethyl-N'-[(8 $\alpha$ )-6-methylergolin-8-yl]-urea  
**Synonyms:** *trans*-dihydro Lisuride, TDHL  
**MF:** C<sub>20</sub>H<sub>28</sub>N<sub>4</sub>O  
**FW:** 340.5  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Terguride is supplied as a solid. A stock solution may be made by dissolving the terguride in the solvent of choice, which should be purged with an inert gas. Terguride is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

Terguride is a receptor modulator.<sup>1</sup> It is an agonist of dopamine D<sub>2</sub> and D<sub>3</sub> receptors and the serotonin (5-HT) receptor subtypes 5-HT<sub>1A</sub>, 5-HT<sub>1B</sub>, and 5-HT<sub>1D</sub>.<sup>2,3</sup> Terguride induces 35SGTPyS binding in CHO cells expressing dopamine D<sub>2</sub> or D<sub>3</sub> receptors (EC<sub>50</sub>s = 0.46 and 0.66 nM, respectively, for the human receptors) or expressing 5-HT<sub>1A</sub>, 5-HT<sub>1B</sub>, or 5-HT<sub>1D</sub> (EC<sub>50</sub>s = 57.54, 426.58, and 14.13 nM, respectively, for the human receptors).<sup>2,3</sup> It is also an antagonist of human  $\alpha$ -adrenergic receptors ( $\alpha$ -ARs; K<sub>i</sub>s = 3.55, 34.67, 3.89, 0.3, 0.45, and 0.76 nM for  $\alpha_{1A}$ ,  $\alpha_{1B}$ ,  $\alpha_{1D}$ ,  $\alpha_{2A}$ ,  $\alpha_{2B}$ , and  $\alpha_{2C}$ -ARs, respectively),  $\beta_1$ - and  $\beta_2$ -ARs (K<sub>i</sub>s = 660.7 and 19.95 nM, respectively), the histamine H<sub>1</sub> receptor (K<sub>i</sub> = 338.84 nM), 5-HT<sub>2D</sub>, 5-HT<sub>2A</sub>, 5-HT<sub>2B</sub>, and 5-HT<sub>2C</sub> receptors (K<sub>i</sub>s = 16.22, 4.79, 7.08, and 47.86 nM, respectively), and dopamine D<sub>1</sub>, D<sub>4</sub>, and D<sub>5</sub> receptors (K<sub>i</sub>s = 28.18, 8.13, and 7.63 nM respectively). Terguride inhibits 5-HT-induced vasoconstriction in rat isolated and perfused lungs when used at concentrations of 1, 3, or 10  $\mu$ M.<sup>4</sup> It inhibits monocrotaline-induced increases in pulmonary blood pressure, cardiac hypertrophy, and decreases in oxygen gas exchange in a rat model of pulmonary hypertension when administered at doses of 0.4 or 1.2 mg/kg per day. Formulations containing terguride have been used in the treatment of hyperprolactinemia.

### References

1. Millan, M.J., Maiofiss, L., Cussac, D., *et al.* Differential actions of antiparkinson agents at multiple classes of monoaminergic receptor. I. A multivariate analysis of the binding profiles of 14 drugs at 21 native and cloned human receptor subtypes. *J. Pharmacol. Exp. Ther.* **303**(2), 791-804 (2002).
2. Newman-Tancredi, A., Cussac, D., Audinot, V., *et al.* Differential actions of antiparkinson agents at multiple classes of monoaminergic receptor. II. Agonist and antagonist properties at subtypes of dopamine D<sub>2</sub>-like receptor and  $\alpha_1/\alpha_2$ -adrenoceptor. *J. Pharmacol. Exp. Ther.* **303**(2), 805-814 (2002).
3. Newman-Tancredi, A., Cussac, D., Quentric, Y., *et al.* Differential actions of antiparkinson agents at multiple classes of monoaminergic receptor. III. Agonist and antagonist properties at serotonin, 5-HT<sub>1</sub> and 5-HT<sub>2</sub>, receptor subtypes. *J. Pharmacol. Exp. Ther.* **303**(2), 815-822 (2002).
4. Dumitrascu, R., Kulcke, C., Königshoff, M., *et al.* Terguride ameliorates monocrotaline-induced pulmonary hypertension in rats. *Eur. Respir. J.* **37**(5), 1104-1118 (2011).

#### 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, 11/04/2024

#### 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