

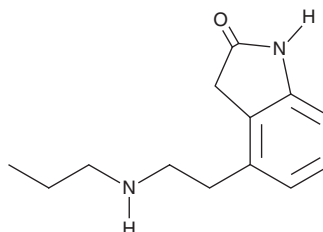
# PRODUCT INFORMATION



## N-Despropyl Ropinirole

Item No. 34156

**CAS Registry No.:** 106916-16-9  
**Formal Name:** 1,3-dihydro-4-[2-(propylamino)ethyl]-2H-indol-2-one  
**Synonym:** SK 104557  
**MF:** C<sub>13</sub>H<sub>18</sub>N<sub>2</sub>O  
**FW:** 218.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 251 nm  
**Supplied as:** A 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

N-Despropyl ropinirole is supplied as a solid. A stock solution may be made by dissolving the N-despropyl ropinirole in the solvent of choice, which should be purged with an inert gas. N-Despropyl ropinirole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of N-despropyl ropinirole in these solvents is approximately 1, 30, and 25 mg/ml, respectively.

N-Despropyl ropinirole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-despropyl ropinirole should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. N-Despropyl ropinirole has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

N-Despropyl ropinirole is an active metabolite of the dopamine D<sub>2</sub> receptor agonist ropinirole (Item No. 23871).<sup>1</sup> It increases extracellular acidification rates in CHO cells expressing human dopamine D<sub>2</sub>, D<sub>3</sub>, and D<sub>4</sub> receptors (EC<sub>50</sub>s = 0.63, 0.063, and 1.23 μM, respectively). N-Despropyl ropinirole is also a potential impurity found in commercial preparations of ropinirole.<sup>2</sup>

### References

1. Coldwell, M.C., Boyfield, I., Brown, T., *et al.* Comparison of the functional potencies of ropinirole and other dopamine receptor agonists at human D<sub>2(long)</sub>, D<sub>3</sub> and D<sub>4.4</sub> receptors expressed in Chinese hamster ovary cells. *Br. J. Pharmacol.* **127**(7), 1696-1702 (1999).
2. Coufal, P., Stulik, K., Claessens, H.A., *et al.* Separation and quantification of ropinirole and some impurities using capillary liquid chromatography. *J. Chromatogr. B Biomed. Sci. Appl.* **732**(2), 437-444 (1999).

#### 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, 11/01/2022

#### 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](http://WWW.CAYMANCHEM.COM)