

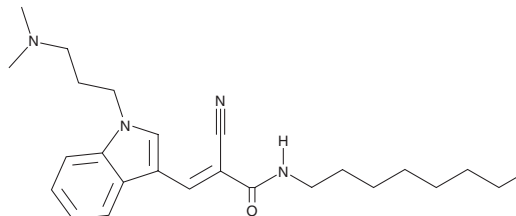
# PRODUCT INFORMATION



## Dynole 34-2

Item No. 34073

**CAS Registry No.:** 1128165-88-7  
**Formal Name:** 2-cyano-3-[1-[3-(dimethylamino)propyl]-1H-indol-3-yl]-N-octyl-2-propenamide  
**MF:** C<sub>25</sub>H<sub>36</sub>N<sub>4</sub>O  
**FW:** 408.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 274, 381 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

Dynole 34-2 is supplied as a solid. A stock solution may be made by dissolving the dynole 34-2 in the solvent of choice, which should be purged with an inert gas. Dynole 34-2 is soluble in organic solvents such as ethanol and DMSO. The solubility of dynole 34-2 in these solvents is approximately 50 and 100 mM, respectively.

### Description

Dynole 34-2 is an inhibitor of dynamin (IC<sub>50</sub>s = 6.9 and 14.2 μM for dynamin 1 and dynamin 2 GTPase activity, respectively).<sup>1</sup> It induces multinucleation, inhibits cytokinesis, and induces apoptosis in HeLa cells when used at a concentration of 10 μM. Dynole 34-2 inhibits the proliferation of several cancer cell lines (GI<sub>50</sub>s = 5.8-15 μM).

### Reference

1. Chircop, M., Pera, S., Mariana, A., *et al.* Inhibition of dynamin by dynole 34-2 induces cell death following cytokinesis failure in cancer cells. *Mol. Cancer Ther.* **10(9)**, 1553-1562 (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, 01/03/2023

#### 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