

# ML60218

## RNA Polymerase III inhibitor

A broad spectrum inhibitor of RNA polymerase III from *S. cerevisiae* as well as from human cells. Combined treatment with ML60218 and an HDAC inhibitor caused augmented suppression of colony formation, induction of cell cycle arrest and apoptotic cell death in human pancreatic adenocarcinoma cells.

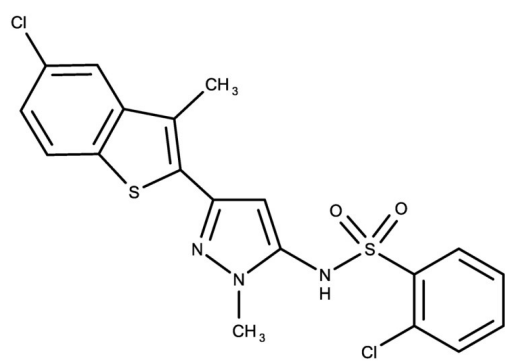
### Ordering Information

[Order Online »](#)

|                 |      |
|-----------------|------|
| ENZ-CHM485-0025 | 25mg |
|-----------------|------|

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** Off-white solid.

**CAS** 577784-91-9

**Couple Target** RNA polymerase

**Couple Type** Inhibitor

**Formula**  $C_{19}H_{15}Cl_2N_3O_2S_2$

**Identity** Determined by NMR.

**MW** 452.38

**Purity** ≥98% (TLC)

**Solubility** Soluble in DMSO (35 mg/ml) ethanol (5 mg/ml).

**Last modified: February 27, 2025**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)