

# XAV939

## PARP5a inhibitor

Potent inhibitor of Tankyrase 1 ( $IC_{50}=11nM$ ) and Tankyrase 2 ( $IC_{50}=4nM$ ). Inhibition of tankyrase promotes stabilization of axin resulting in  $\beta$ -catenin degradation and Wnt signalling disruption. Inhibits proliferation of the  $\mu$ -catenin-dependent colon carcinoma cell line DLD-1. Promotes cardiomyogenic development in mesoderm progenitor cells.

Citations: 3

[View Online »](#)

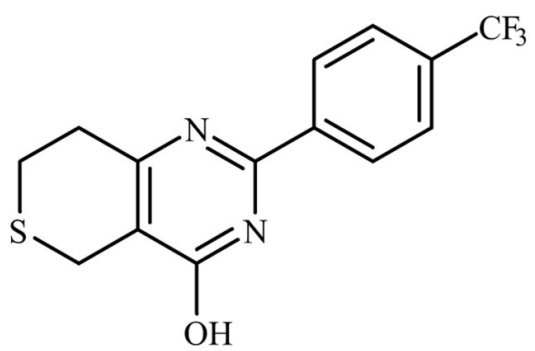
## Ordering Information

[Order Online »](#)

|                |      |
|----------------|------|
| BML-WN100-0005 | 5mg  |
| BML-WN100-0025 | 25mg |

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |                                                                  |
|-------------------|------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. |
| Handling          | Solutions can be stored at -20°C for up to 3 months.             |
| Long Term Storage | Ambient                                                          |
| Shipping          | Ambient Temperature                                              |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                                      |
|------------------|--------------------------------------------------------------------------------------|
| Alternative Name | 3,5,7,8-Tetrahydro-2-[4-(trifluoromethyl)phenyl]-4H-thiopyrano[4,3-d]pyrimidin-4-one |
| Appearance       | White solid.                                                                         |
| CAS              | 284028-89-3                                                                          |
| Couple Target    | PARP                                                                                 |
| Couple Type      | Inhibitor                                                                            |
| Formula          | $C_{14}H_{11}F_3N_2OS$                                                               |
| MW               | 312.3                                                                                |
| Purity           | ≥98% (HPLC)                                                                          |
| Solubility       | Soluble in DMSO (20 mg/ml).                                                          |

Last modified: May 29, 2024



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-fr@enzolifesciences.com](mailto:info-fr@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)