

## ML323

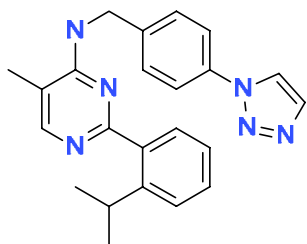
Cat. # SI9634

**Background:** ML323 binds to the complex of USP1-UAF1 resulting in the inhibition of USP1 deubiquitylase activity (Liang, Dexheimer et al. 2014). The  $IC_{50}$  is reported to be 76nM against Ub-Rho110 and 174nM against K63-linked diubiquitin. Inhibition is reversible and exhibits mixed-mode kinetics. ML323 showed little or no inhibitory activity against a collection of DUBs including USPs 2, 5, 7, 8, 10, 11, 14, and 21, UCH-L1 and UCH-L3, and SenP1.

**Application:** Studying the role of USP1-UAF1 in DNA-damage repair. Sensitization of cells to cisplatin-killing.

### Product Information

<b>CAS No.:</b>	1572414-83-5
<b>Purity:</b>	99%, NMR consistent with structure
<b>Molecular Weight:</b>	384.48
<b>Physical State:</b>	Powder
<b>Quantity:</b>	5mg
<b>Solubility:</b>	198mM in DMSO
<b>Storage:</b>	Store desiccated as supplied at -20°C for 2 years
<b>Formula:</b>	$C_{23}H_{24}N_6$
<b>SMILES String:</b>	<chem>Cc1cnc(nc1NCc2ccc(cc2)n3ccnn3)c4cccc4C(C)C</chem>



### References

Liang, Q., T. S. Dexheimer, et al. (2014). "A selective USP1-UAF1 inhibitor links deubiquitination to DNA damage responses." *Nat Chem Biol* **10**(4): 298-304.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2013 LifeSensors, Inc. All Rights Reserved