

SKPin C1

2-[4-Bromo-2-[[4-oxo-3-(3-pyridinylmethyl)-2-thioxo-5-thiazolidinylidene]methyl]phenoxy]acetic acid

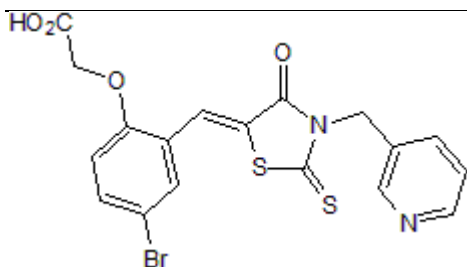
Cat. # SI9845

Background: Inhibits the cullin-RING ubiquitin E3 ligase SCF-Skp2. Selectively inhibits Skp2-mediated p27 degradation by reducing p27 binding. In cancer cells the compound induces p27 accumulation and promotes cell type-specific blocks in G1 or G2/M phases^{1,2}.

Application: Cell-based and *in vitro* studies of Skp2 function

Product Information:

CAS No.	432001-69-9
Purity:	98% (TLC)
Molecular Weight:	465.34 Da
Physical State:	Powder
Quantity:	5mg, 25mg
Solubility:	DMSO (58mg/ml); Ethanol (8mg/ml); Water (<1mg/ml)
Storage:	Store desiccated as supplied at 4°C.



Formula:

C₁₈H₁₃BrN₂O₄S₂

References

1. Rico-Bautista, E. and D. A. Wolf (2012). "Skipping cancer: small molecule inhibitors of SKP2-mediated p27 degradation." *Chem Biol* **19**(12): 1497-1498.
2. Wu, L., A. V. Grigoryan, et al. (2012). "Specific small molecule inhibitors of Skp2-mediated p27 degradation." *Chem Biol* **19**(12): 1515-1524.

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