

SL-01

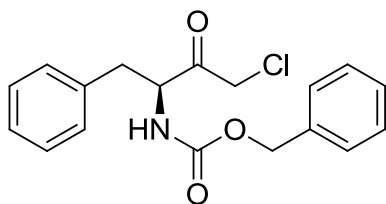
N-Benzyloxycarbonyl-L-phenylalaninylchloromethyl ketone; ZPCK

Cat. # SI9860

Background: Inhibits the p53-MDM2 interaction (20 μ M). Inhibits U2Os cell growth (EC₅₀=10 μ M). Has also been shown to inhibit bovine chymotrypsin A- γ .

Product Information:

CAS No.	26049-94-5
Purity:	>98% (TLC); NMR (Conforms)
Molecular Weight:	331.8
Physical State:	White Crystalline Solid
Quantity:	50 mg
Solubility:	May be dissolved in DMSO (50 mg/ml) or Ethanol (15mg/ml, warm)
Storage:	Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.



Formula: C₁₈H₁₈ClNO₃

References

1. Li J. *et al.* A cell-based high-throughput assay for the screening of small-molecule inhibitors of p53-MDM2 interaction. *J. Biomol. Screening* 2011 **16**:450

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2011 LifeSensors, Inc. All Rights Reserved