

## (-)-Blebbistatin \*CAS#: 856925-71-8\*

Catalog number: 80001 Unit size: 1 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	24 months upon receiving
Chemical Properties	
Appearance	Yellow solid
Molecular Weight	292.33
Soluble In	DMSO
Chemical Structure	

## Applications

(-)-Blebbistatin is a cell-permeable inhibitor for non muscle myosin II ATPase with IC50 of 2  $\mu$ M. It does not inhibit myosin light chain kinase, inhibits contraction of the cleavage furrow without disrupting mitosis or contractile ring assembly.