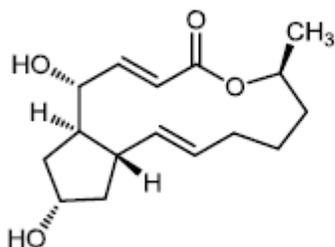


Brefeldin A

Cat. # AC-2007

Background: Brefeldin A is a fermentation product from *Penicillium brefeldianum*. It is a lactone antibiotic and ATPase inhibitor for protein transport with IC_{50} of 0.2 μ M in HCT 116 cells. Brefeldin A has been widely used as specific inhibitor of protein translocation from endoplasmic reticulum to Golgi¹. Blocks secretion and prevents the assembly of non-clathrin-coated buds on Golgi cisternae². Inhibits guanine nucleotide-exchange proteins BIG1 and BIG2, activators of ADP-ribosylation factors³. Induces apoptosis independently of p53⁴. Activates sphingomyelin metabolism⁵. Cell permeable.

Application: Autophagy inducer



Product Information

Purity:	>99.0% by HPLC
MW:	280.4
Formula:	C ₁₆ H ₂₄ O ₄
CAS No.	20350-15-6
Physical State:	Lyophilized powder
Amount:	5mg, 25mg, 50mg
Solubility:	100 mg/mL in DMSO; or 25 mg/ml in warm ethanol
Storage:	Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.

References

1. T Fujiwara *et al.* *J. Biol. Chem.* 1988 263:18545
2. L Orci *et al.* *Cell* 1991 64:1183
3. KF Xu *et al.* *Proc. Natl. Acad. Sci. USA* 2005 102:2784
4. RG Shao *et al.* *Exp. Cell Res.* 1996 227:190
5. CM Linardic *et al.* *Cell Growth Differ.* 1996 7:765

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