

PRODUCT DATA SHEET



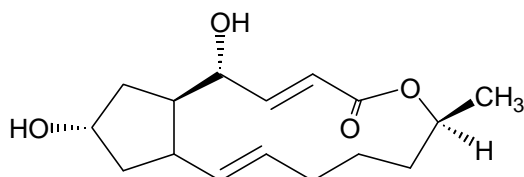
PRODUCT: Brefeldin A
BFA

CAS NO: 20350-15-6

CATALOG NO.: G-405

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 204-205°C)

MOLECULAR FORMULA: C₁₆H₂₄O₄

MOLECULAR WEIGHT: 280.4

PURITY: 98%(TLC)

SOLUBILITY: Soluble in ethanol (5 mg/mL) and DMSO (20 mg/mL)

STORAGE: Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

APPLICATION NOTES: Inhibits binding of the cytosolic coat protein, β-COP and ARF to Golgi membranes.¹⁻³ Induces a rapid redistribution of Golgi into the ER and blocks transport of proteins into post-Golgi compartments.⁴ Blocks the action of ARF activating protein.⁵

- REFERENCES:
1. J.G.Donaldson *et al.* *Science* 1991 **254** 1197
 2. J.G.Donaldson *et al.* *J.Cell Biol.* 1991 **112** 579
 3. L.Orci *et al.* *Cell* 1991 **64** 1183
 4. L.Lippincott-Schwartz *et al.* *Cell* 1990 **60** 821
 5. P.A.Randazzo *et al.* *J.Biol.Chem.* 1993 **268** 9555
 6. J.G.Donaldson *et al.* *Nature* 1992 **360** 350
 7. J.B.Helms *et al.* *Nature* 1992 **360** 352
 8. P. Chardin and F. McCormick *Cell* 1999 **97** 153

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.