

Okadaic acid, free acid

Phosphatase Inhibitor

Catalog # P16-902

Lot # P3210-1

CAS #78111-17-8

Product Description

Molecular Formula: C₄₄H₆₈O₁₃

Appearance: White lyophilized solid

Melting Point: N/A

Molecular Weight: 805

Purity: >98% (TLC)

Solubilization: May be dissolved in DMSO (40 mg/ml); or Ethanol (5 mg/ml)

Alias

Specific Activity

PP2A and PP1 inhibitor

Storage and Stability

Store desiccated as supplied at -20°C for up to 2 years.

Store solutions at -20°C for up to 1 month.

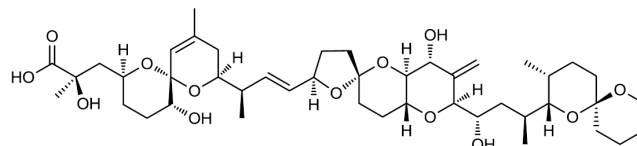
Scientific Background

A naturally occurring polyether marine toxin. Potent and selective inhibitor of protein phosphatases inhibiting PP2A completely at 1nM and PP1 at higher concentrations, while PP2C is not inhibited. It is an extremely useful tool for studying cellular protein phosphorylation (2). Induces apoptosis in a variety of cell types (3). Activates atypical protein kinase isotypes and induces insulin-like effects in adipocytes (4).

References

1. P Cohen *et al.* *Trends Biochem. Sci.* 1990 15:98.
2. DG Hardie *et al.* *Methods Ezymol.* 1991 201:469.
3. TJ Haneji *et al.* *Int. J. Oncol.* 2013 42:1904.
4. ML Standaert *et al.* *J. Biol. Chem.* 1999 274:14074.

Molecular Structure



Okadaic acid, free acid

Phosphatase Inhibitor

Catalog # P16-902

Lot # P3210-1

CAS # 78111-17-8

Purity >98% (TLC)

Format Lyophilized powder

Solubilization May be dissolved in DMSO (40 mg/ml); or Ethanol (5 mg/ml).

Storage & Shipping Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.