

Okadaic Acid

Catalog #: 27047

Lot #: 181008

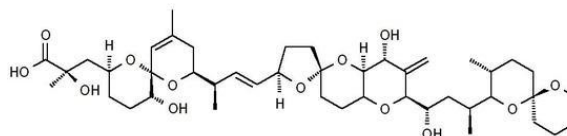
Size: 100 µg

CAS Registry #: 78111-17-8

Structure:

Purity: ≥98%

Chemical Formula: C₄₄H₆₈O₁₃



Molecular Weight: 805.0

Description: Potent inhibitor of protein phosphatases, especially the PP-1 and PP-2 classes, in numerous cell types. Non-phorbol type tumor promoter.

Synonym(s): Halochondrine A

Appearance: Off-white solid.

Solubility: Soluble in DMSO or ethanol.

Biological Activity: Okadaic acid increases intracellular protein phosphorylation and acts as a potent tumor promoter.

Storage/Stability: Store at or below -20°C. Solid form is stable at least 12 months from date of receipt, when stored as directed. Do not store aqueous solutions for more than one day.

Warnings: Highly irritant to skin and mucous membranes. Tumor promoter. Gloves and mask should be worn when using this compound. Care must be taken to prevent contact through all routes of exposure.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Cohen, P. *et al.*, *Trends Biochem. Sci.* 1990;**15**, 98-102.
2. Saganuma, M. *et al.*, *Proc. Nat. Sci. USA.* 1988;**85**, 1768-1771.