

---

## Ginkgolic acid

(Catalog No. M75004)

---

### Alternative Names

(Z)-6-(Pentadec-8-enyl)-2-hydroxybenzoic acid

### Description

Cell-permeable anacardic acid analog that inhibits protein SUMO modification (IC<sub>50</sub> = 3.0 μM using RanGAP1-C2 as substrate) in an ATP-dependent manner by selectively targeting SUMO-activating enzyme E1 and interfering with E1-SUMO intermediate formation independent of E1 active site cysteine.

### CAS Number

22910-60-7

### Chemical Formula

C<sub>22</sub>H<sub>34</sub>O<sub>3</sub>

### Molecular Weight

346.5

### Size

5 mg

### Purity

≥80%

### Solubility

DMSO or EtOH

### Storage Temp

-20°C

### Ordering Information:

Products	Size	Cat. No.
Ginkgolic acid	5 mg	M75004 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*