

## Z-VAD-FMK

**Catalog #:** 27313

**Lot #:** 170620

**Size:** 5 mg

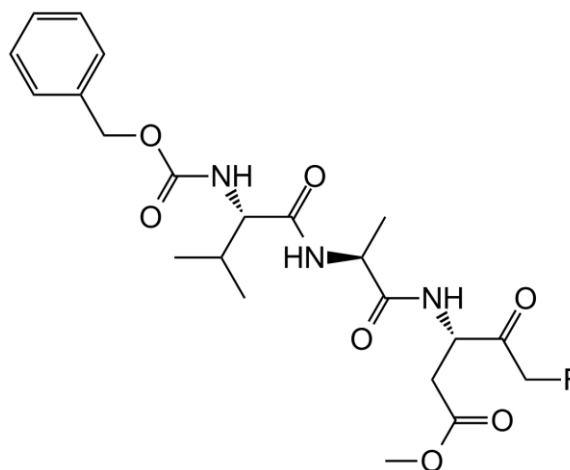
**CAS Registry #:** 187389-52-2

**Structure:**

**Purity:** ≥97%

**Chemical Formula:** C<sub>22</sub>H<sub>30</sub>FN<sub>3</sub>O<sub>7</sub>

**Molecular Weight:** 467.49



**Description:** Z-VAD(OMe)-FMK is a cell permeable, irreversible pan-caspase inhibitor and anti-apoptotic agent. It has been shown to inhibit Fas-mediated apoptosis in Jurkat T cells and apoptosis in THP.1 cells induced by diverse stimuli (including cycloheximide, thapsigargin, etoposide and staurosporine). Inhibits caspases 1, 3, 4, and 7.

**Storage:** Store at -20°C

**Solubility:** DMSO (>10 mg/ml), dimethyl formamide (7 mg/ml) or 100% ethanol (1.3 mg/ml)

### References:

1. Rodríguez-Enfedaque, A., et al. *Biochim. Biophys. Acta.* 1823: 1343-1352 (2012).