

Z-FA-FMK: Caspase Inhibitor Negative Control

Catalog No.: NBP2-29384, 3 mg

NBP2-29384, 5 mg

Sequence: Z-Phe-Ala-FMK
Z-FA-FMK

Molecular Weight: 386

Storage: Key cool and dry. The solid product is stable in the dessicator at room temperature for 1 year. However, we recommend storing dessicated at -20°C.

Form: White Solid

Analytical Data:

Mass Spec: M+1=387.0

Chromatography: TLC:EtOAc: 4, Hex: 6, Single Spot, Rf:0.4

H NMR: All functional groups are present

Solubility: Make a stock solution of 10 mM in high purity DMSO (>99.9%). The stock solution is stable at -20°C for 6-8 months. Avoid repeated freeze/thaw cycles of the stock solution. For multiple uses, we suggest aliquoting the stock solution prior to freezing. Bring the solution to room temperature before opening the vial cap.

Description:

Z-FA-FMK is an inhibitor of cathepsins B and L. Z-FA-FMK is designed for *in vitro* use as a cell permeable control for the effects of fluormethyl-ketone derivatives (FMKs) on caspase activity; it has been used as a negative control for FMK based inhibitors of caspases in a various studies (1-4).

References:

1. Sillence, D.J. and D. Allan. 1997. Biochem. J. 324:29
2. Gottron, F.J. et al. 1997. Mol. Cell. Neurosci. 9:159.
3. McColl, K.S. et al. 1998. Mol. Cell. Endocrinol. 139:229
4. Gregoli, P.A. and M.C. Bondurant. 1999. J. Cell Physiol. 178:133.

Assay Method for using Z-FA-FMK as a Negative Control for Caspase Inhibitors

Materials:

Dissolve 1mg of in 259 μ l DMSO for a 10 mM solution.

Method:

Add 2 μ l of 10 mM stock solution in 1 ml of culture medium = 20 μ M final Z-FA-FMK concentration and final DMSO concentration of 0.2%. Levels of DMSO above this may cause some cellular toxicity thus masking the effect of the ICE-protease inhibitors.

Caspase FMK inhibitors such as Z-VAD-FMK, BOC-D-FMK, Z-DEVD-FMK, Z-IETD-FMK, and Z-LEHD-FMK are generally used in tissue culture to inhibit apoptosis at final working concentrations of 50 nM -100 μ M. Optimal concentrations may vary due to cell type, method of apoptosis induction, and length of culture. Each investigator should determine the most effective inhibitor concentration for their particular assay. The Z-FA-FMK negative control inhibitor should then be used at the same concentration as the caspase inhibitor.