



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP2-29396

NBP2-29396 Protocol

Product Handling - Assay Method (NBP2-29396)

Product Handling - Assay Method (NBP2-29396): https://www.novusbio.com/products/z-d-ome-e-ome-vd-ome-fmk-inhibitor_nbp2-29396

Materials:

Dissolve 1mg of Z-D(OMe)E(OMe)VD(OMe)-FMK in DMSO to get appropriate concentration:

75 ul DMSO = 20mM

150 ul DMSO = 10 mM

300 ul DMSO = 5 mM

Method:

Add 2 ul of above stock solutions to 1 ml of culture medium containing cells to give 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. Two microliter of 10mM stock solution in 1 ml medium = 20 uM final Z-D(OMe)E(OMe)VD(OMe)-FMK concentration. Caspase inhibitors have been successfully used in tissue cultures to inhibit apoptosis at final working concentrations of approximately 50 nM to 100 uM. The variability reflects cell type and method. Investigators should determine the optimal concentration empirically.