



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP2-29397

NBP2-29397 Protocol

Product Handling - Assay Method (NBP2-29397)

Product Handling - Assay Method (NBP2-29397): https://www.novusbio.com/products/z-ie-ome-td-ome-fmk-inhibitor_nbp2-29397

Materials:

Dissolve 5 mg of Z-IETD-FMK in DMSO to get appropriate concentration:

765 ml DMSO = 10 mM

383 ml DMSO = 20 mM

1530 ml DMSO = 5 mM

Method:

Add 2 ml 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. 2 ml of 10 mM stock in 1 ml medium = 20 mM final Z-IETD-FMK concentration.