



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP2-29398

NBP2-29398 Protocol

Product Handling - Assay Method (NBP2-29398)

Product Handling - Assay Method (NBP2-29398): https://www.novusbio.com/products/z-le-ome-hd-ome-fmk-inhibitor_nbp2-29398

Materials:

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

310 ul DMSO = 20 mM

620 ul DMSO = 10 mM

1240 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. 2 ul of 10 mM stock in 1 ml medium = 20 uM final Z-LEHD-FMK concentration.