



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP2-29392

NBP2-29392 Protocol

Product Handling - Assay Method (NBP2-29392)

Product Handling - Assay Method (NBP2-29392): https://www.novusbio.com/products/z-val-ala-asp-ome-fmk-inhibitor_nbp2-29392

Materials:

Dissolve 1mg of Z-Val-Ala-Asp(OMe)-FMK in DMSO to get appropriate concentration:

107 ul DMSO = 20mM

214 ul DMSO = 10 mM

428 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.