



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP2-29395

NBP2-29395 Protocol

Product Handling - Assay Method (NBP2-29395)

Product Handling - Assay Method (NBP2-29395): https://www.novusbio.com/products/boc-d-ome-fmk-inhibitor_nbp2-29395

Materials:

Dissolve 1mg of BOC-D-FMK in DMSO to get appropriate concentration:

190 μ l DMSO = 20mM

381 μ l DMSO = 10 mM

762 μ l DMSO = 5 mM

Method:

Add 2 μ l 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 μ l of 10 mM stock in 1 ml medium = 20 μ M final BOC-D-FMK concentration.