



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

TCM-ProteaseArrest™

A Complete Protease Inhibitor
for Tissue Culture Media

(Cat. # 786-238, 786-239)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

TCM-ProteaseArrest™ is a mixture of protease inhibitors with a broad specificity for the inhibition of serine, cysteine, aspartic (acid) proteases and aminopeptidases. The protease inhibitor cocktail has been developed to be used specifically as a supplement for tissue culture medium to prevent the degradation of secreted proteins.

TCM-ProteaseArrest™ is supplied as a 200X concentrated solution in sterile filtered DMSO. It is highly recommended to determine the appropriate dilution factor for a specific cell line, starting at a 1:200 dilution of TCM-ProteaseArrest™ to prevent harmful effects of DMSO to cell growth.

ITEM(S) SUPPLIED

Description	Cat. # 786-238	Cat. # 786-239
TCM-ProteaseArrest™ [200X]	1ml	2ml

STORAGE CONDITIONS

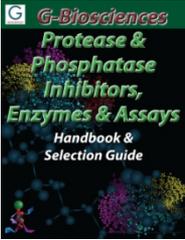
Shipped at ambient temperature, upon arrival, store it at -20°C, and is stable for up to 1 year.

PROTOCOL

1. Allow the solution to warm to room temperature. The solution is in suspension form, vortex or tap the vial before removing the solution.
2. It is recommended to determine the dilution of the protease inhibitor cocktail for each cell line. Start the dilutions at 1:200 as high concentrations may result in a deleterious effect by the DMSO. Further dilution of the cocktail (1:400, 1:800) may be necessary as different cell lines will differ in their sensitivity to the protease inhibitors.
3. The cocktail will remain effective for 24-48 hours depending on the cell line and the culture medium. After this period the media should be replaced with freshly prepared media supplemented with TCM-ProteaseArrest™.

RELATED PRODUCTS

Download our Protease & Phosphatase Inhibitors, Enzyme & Assays Handbook.



<http://info.gbiosciences.com/protease-phosphatase-inhibitors-enzymes-assay-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 05/5/2014 CMH/IA



www.GBiosciences.com