

Protease Inhibitor Cocktail**Cat. # PIC02****Lot# 123****Amount when reconstituted: 1ml of 100X stock****Store at 4°C desiccated and at 4°C or -70°C when reconstituted****Material**

Protease Inhibitor Cocktail is supplied as a white lyophilized powder. The cocktail contains the following protease inhibitors, Pepstatin A, Leupeptin, Benzamidine, Na-p-tosyl-L-arginine methyl ester (TAME). At 1X final working concentration the inhibitors are at 1 µM, 1.5 µM, 1 mM and 0.4 mM respectively.

Mode of Action

pepstatin A	Reversibly inhibits some aspartic proteases
Leupeptin	Reversibly inhibits trypsin like serine and some cysteine proteases
Benzamidine	Inhibits serine proteases
TAME	Inhibits arginine proteases, a thrombin and trypsin inhibitor

Storage and Reconstitution

Lyophilized Protease Inhibitor Cocktail is stable at 4°C desiccated (<10% humidity) for 1 year. Each tube of Protease Inhibitor Cocktail should be resuspended in 1ml of DMSO (dimethyl sulfoxide) to give a 100X strength stock. The reconstituted stock can be stored at 4°C for up to 6 months or at -70°C for up to 1 year.

Uses

Used at 1X final concentration as an inhibitor of a wide range of proteases.

Related Products

See www.cytoskeleton.com for a complete listing.