

PRODUCT NAME		
Protease Inhibitor Mix (P.I.)		
ChIP grade reagent Separately available		
Cat. No: C12010011	Size: 100 µl	Concentration: 200 x stock solution
Cat. No: C12010012	Size: 300 µl	Concentration: 200 x stock solution

Description: This protease inhibitor Mix (P.I.) has been extensively validated in chromatin immunoprecipitation assays (ChIP). It contains inhibitors with a broad specificity for serine, cysteine, and acid proteases, and aminopeptidase.

Our protease inhibitor Mix is used during the shearing and the immunoprecipitation of the chromatin (ChIP). It is included in our OneDay ChIP KIT (cat# kch-onedIP-060, -180). It is now separately available. For more details about these kits, see instruction manual on our web site (www.diagenode.com).

Format: Supplied as a solution.

Storage: For long storage, aliquot and store at -20°C. The product is stable for 1-2 years when stored at -20 °C and 4 months at 4°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Availability date: October 29, 2007

Last data sheet update: October 26, 2007