

PRODUCT NAME			
Protease Inhibitor Mix (P.I.)			
ChIP grade reagent Separately available			
Cat. No: <b>kch-107-020</b>	Size: 20 µl	Concentration: 200 x stock solution	Lot #: PI103

**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](http://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

**Lot #:** PI103/ aliquoting day: October 02, 2007