

| PRODUCT NAME | | | |
|--|----------------|---------------------|-------------|
| Protease Inhibitor Mix (P.I.) | | | |
| ChIP grade reagent Separately available | | | |
| Cat. No: kch-405-001 | Size: 1 tablet | Concentration: 25 x | Lot #: D002 |

Description: This protease inhibitor Mix (P.I.) has been extensively validated in chromatin immunoprecipitation assays (ChIP). It inhibits a broad spectrum of serine, cysteine and metalloproteases as well as calpains.

Our protease inhibitor Mix is used during the immunoprecipitation of the chromatin (ChIP). It is included in our **Transcription ChIP KIT (cat# kch-redTBP-012)** and our **Histone ChIP KIT (cat# kch-orgHIS-012)**. 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 tablet. Dissolve the tablet of protease inhibitor mix (P.I.) in 400 µl of water. It is then a 25x stock solution.

Storage: The tablet is stable at -20°C for at least 2 years. The stock solution is stable at -20°C for at least 4 months. 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 #: D002/ aliquoting day: October 02, 2007