

Protease Inhibitor Mix (P.I.)

ChIP grade reagent
Separately available

Cat. No. **C12010015** Format: 1 tablet

Last datasheet update: August 29, 2019

Product 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.

The protease inhibitor mix is recommended during chromatin shearing and ChIP incubation.

Format

Supplied as a tablet. Dissolve the tablet of protease inhibitor mix (P.I.) in 400 μ l of water to obtain a 25x stock solution.

Storage and stability

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.