

## Protease Inhibitor Cocktail

---

**Catalog No:** 37490, 37491

**Format:** 1 ml, 5 ml

### Quality Control:

Active Motif's Protease Inhibitor Cocktail is a broad-spectrum cocktail designed to prevent proteolytic degradation during lysis and extraction procedures from mammalian cells or tissues. The cocktail solution inhibits aminopeptidases and cysteine, serine and aspartic acid proteases. Protease Inhibitor Cocktail is provided EDTA-free, as a ready-to-use 100X stock solution in DMSO. The cocktail composition and inhibitor targets are listed below:

- AEBSF, 104 mM - serine proteases
- Aprotinin, 80 µM - serine protease
- Bestatin, 4 mM - aminopeptidases
- E-64, 1.4 mM - cysteine proteases
- Leupeptin, 2 mM - serine and cysteine proteases
- Pepstatin A, 1.5 mM - aspartic acid proteases

It is recommended to use 10 µl of the cocktail solution for every 1 ml of cell lysate. Since protease levels may vary among cell and tissue types, it may be necessary to increase the concentration of inhibitors.

**Storage and Guarantee:** The Protease Inhibitor Cocktail is provided in DMSO as a faint yellow solution. The product is shipped at room temperature and should be stored at -20 °C long-term. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.