

## 100 mM PMSF

---

**Catalog No:** 37495

**Format:** 0.5 ml

**Quality Control:** Active Motif's 100 mM PMSF (part number 101664), also known as Phenylmethylsulfonyl fluoride, is a non-specific inhibitor of serine proteases and the cysteine protease papain (reversible by DTT treatment). PMSF does not inhibit metalloproteases, most cysteine proteases or aspartic proteases.

PMSF is provided at a concentration of 100 mM in isopropanol. It is recommended for use at a working concentration of 0.1 - 1 mM. Since protease levels may vary among cell and tissue types, it may be necessary to optimize the concentration of inhibitors.

**Storage and Guarantee:** Store at -20°C. 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.