

PRODUCT DATA SHEET



TIMP-2 (human, purified)

CATALOG NO.: SE-469

LOT NO.: XXXXX

DESCRIPTION: MW approximately 22 kDa by SDS-PAGE. Tissue inhibitor of metalloproteinases-2 (TIMP-2) purified from human fibroblasts. Free of MMPs or other TIMPs.

PURITY: >95% by SDS-PAGE.

USAGE: MMP inhibitor *in vitro*¹ or in tissue culture²; positive control in antibody applications or reverse zymograms. TIMP-2 inhibits most MMPs, particularly MMP-2³⁻⁵. It can also inhibit ADAMs such as ADAMTS⁶. TIMP-2 is also required for activation of pro-MMP-2 by MT1-MMP and MT2-MMP⁷. While it can inhibit angiogenesis, it can also promote cell growth^{7,8}.

SUPPLIED AS: 5 µg at 0.2 µg/µl in 50 mM Tris-HCl pH 7, 200 mM NaCl, 5 mM CaCl₂, 1 µM ZnCl₂, 0.05 % Brij-35, and 0.05% NaAzide.

STORAGE: -80°C. After initial defrost, aliquot product into individual tubes and refreeze at -80°C. Avoid repeated freeze/defrost cycles.

NOTE: When stored under the above conditions, this protein is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing could lead to loss of activity.

REFERENCES:

1. M.W. Olson *et al.* *J. Biol. Chem.* 1997 **272** 29975
2. J. Oh *et al.* *Cancer Res.* 2004 **64** 9062
3. M. Hutton *et al.* *Biochemistry* 1998 **37** 10094
4. H. Nagase and K. Brew *Arthritis Res.* 2002 **4(suppl 3)** S51
5. E. Lambert *et al.* *Crit. Rev. Oncol. Hematol.* 2004 **49** 187
6. G. Hashimoto *et al.* *FEBS Lett.* 2001 **494** 192
7. R. Visse and H. Nagase *Circ. Res.* 2003 **92** 827
8. M. Takigawa *et al.* *Biochem. Biophys. Res. Commun.* 1990 **171** 1264

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.