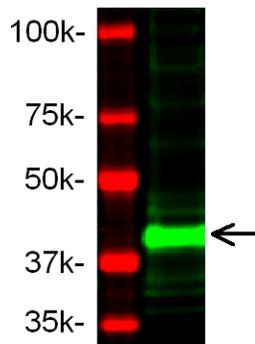


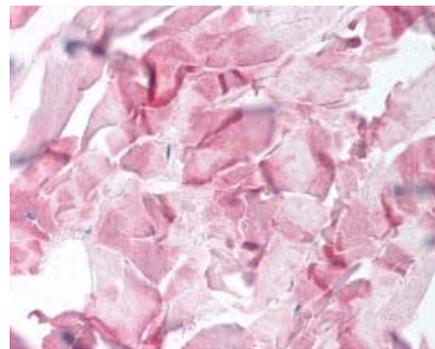


Product Data Sheet

Product Name:	Anti-MMP-1 (NT)
Catalog Number:	29574
Lot Number:	See label on vial
Product Description:	This polyclonal antibody is supplied as peptide affinity purified rabbit antiserum, 50 µg (at 200 µg/ml) in 1X PBS (pH 7.4) containing 0.1% BSA and 0.05% sodium azide.
Immunogen:	Rabbit polyclonal anti-MMP-1 (NT) antibody was raised against a synthetic collagenase-1 peptide derived from the N-terminus of human MMP-1.
Species Reactivity:	This antibody is specific to human.
Application Notes:	The following concentration ranges are recommended starting points for this product. Optimal working concentrations should be determined by the investigator for specific applications.
IHC:	5.0-10.0 µg/ml
WB:	1.0-2.0 µg/ml for 2 hrs at RT



Western blot analysis of MMP-1 in human MMP-1 over-expressing HEK293 cell lysate probed with the anti-MMP-1 (NT) antibody (cat#29574).



Immunocytochemistry of human dermal collagen in the skin (Formalin fixed / Paraffin-embedded) stained with rabbit anti-MMP-1 (NT) (Cat#29574).

Background:

Matrix metalloproteinases (MMP) are proteolytic enzymes capable of degrading connective tissue components. Imbalanced secretion of certain MMP or disturbances in the differential control of MMP by tissue inhibitor of MMPs (TIMPs) is implicated in the invasive potential of malignant tumors. Interstitial collagenase (MMP-1)(52 kDa), gelatinase A/type IV collagenase (MMP-2) (72 kDa), and gelatinase B/type V collagenase (MMP-9) (92 kDa) dissolve extracellular matrix (ECM) and may initiate and promote angiogenesis.

Storage:

Store at 2-8 °C for up to one year. Avoid repeated freezing and thawing.

This product is for *in vitro* research use only.