

Catalog #	Aliquot Size
<b>M846-40N-10</b>	<b>10 µg</b>
<b>M846-40N-100</b>	<b>100 µg</b>
<b>M846-40N-1000</b>	<b>1 mg</b>

## Recombinant Human MMP2

Recombinant human protein expressed in E. coli cells

**Catalog # M846-40N**

Lot # G3211-98

### Product Description

Recombinant Human MMP2 was expressed in E. coli cells. The protein accession number is [P08253](#)

### Gene Aliases

MMP-2, 72 kDa type IV collagenase, 72 kDa gelatinase, Gelatinase A, TBE-1, Recombinant Human Matrix Metalloproteinase-2 (MMP2)

### Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method.

### Formulation

Supplied as a 0.2 µm filtered solution in Tris, CaCl<sub>2</sub>, NaCl and Brij-35

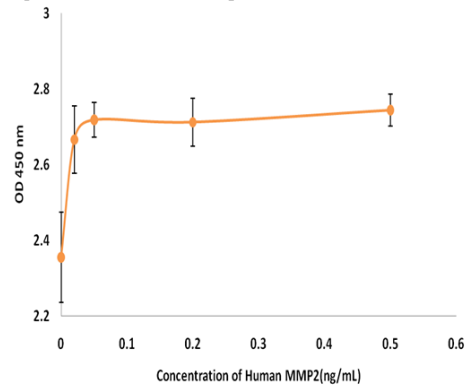
### Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

### Scientific Background

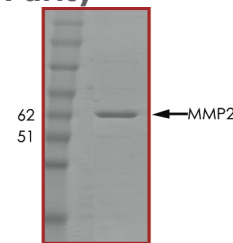
MMP-2 belongs to a family of endoproteases which need calcium and zinc in order to exert their catalytic capability (extracellular matrix remodeling). The elevated expression of these enzymes has been linked to many degenerative disorders, such as arthritis, cancer metastasis and skeletal growth-plate disorders. It has been shown that MMP-2 associates with many connective tissue cells as well as neutrophils, macrophages and monocytes.

### Specific Activity



The ED (50) was determined by the dose-dependent proliferation of immortalized human pulmonary fibroblasts cells and was found to be <0.05ng/mL.

### Purity



Recombinant Human MMP2 resolved on a 10% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **62.0 kDa**

## Recombinant Human MMP2

Recombinant human protein expressed in E. coli cells

Catalog Number	M846-40N
Specific Lot Number	G3211-98
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**