MMP-2 (mouse fibroblasts), (purified)

Ordering Information

Order Online »

BML-SE505-0005

5µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Matrix metalloproteinase 2, Gelatinase A, 72 kDa Type IV

collagenase

Application NotesUseful for the study of enzyme regulation and kinetics,

cleavage of substrates in vitro or in tissue culture,

zymography or immunoblot positive control.

Formulation Liquid. In 50mM Tris-HCl, pH 7.0, containing 200mM NaCl,

5mM CaCl₂, 1µM ZnCl₂, 0.05% sodium azide and 0.05%

Brij 35.

Purity Detail Free of TIMPs and other MMPs.

Source Isolated from mouse fibroblasts. Primarily latent form.

UniProt ID P08253

Last modified: May 29, 2024



eu@enzolifesciences.com