## HGF/SF monoclonal antibody (A10)

HGF/SF (Hepatocyte growth factor/Scatter factor) is a multifunctional heterodimeric polypeptide. It mediates the growth and scattering of various cell types, epithelial mesenchymal transition, the formation of tubules and lumens, and promotes angiogenesis. The ligand-receptor pair has also shown to be uniquely involved in most human solid tumors and to participate in tumor development, invasion, and metastasis.

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

**Ordering Information** 

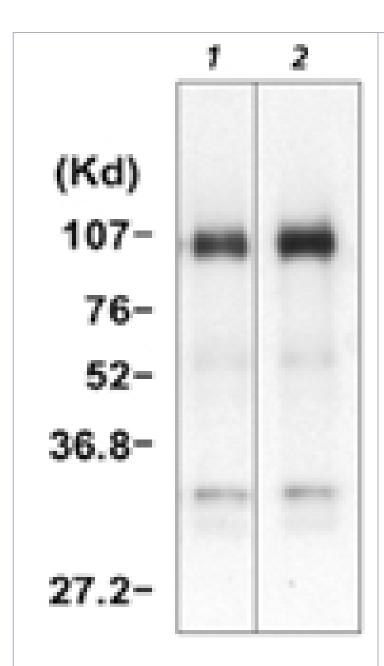
Order Online »

ADI-905-163-100

100µg

Manuals, SDS & CofA

**View Online »** 



Immunoprecipitation analysis: Lane 1 anti-human HGF Clone A10 and Lane 2 positive anti-serum.

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name HGFB, HPTA, hepatocyte growth factor

**Application** IP, WB

Clone A10

**Formulation** Liquid. In PBS containing 0.01% sodium azide.

Concentration 1mg/ml.

**Host** Mouse

Immunogen Human Hepatocyte Growth Factor/Scatter Factor.

lsotype lgG1

Purity Detail Protein G affinity purified.

**Source** Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

UniProt ID P14210

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

