

## Product datasheet

# Mouse HGF ELISA Kit ab100687

5 References 1 Image

### Overview

<b>Product name</b>	Mouse HGF ELISA Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture extracts, Tissue Extracts
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 100 pg/ml
<b>Range</b>	0.25 ng/ml - 60 ng/ml
<b>Recovery</b>	100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture extracts	114.4	100% - 134%
Tissue Extracts	120.2	110% - 130%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Mouse

**Product overview** Abcam's HGF Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse HGF in cell lysates and tissue lysates.

This assay employs an antibody specific for mouse HGF coated on a 96- well plate. Standards and samples are pipetted into the wells and HGF present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse HGF antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of HGF bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Notes** Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

**Platform** Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

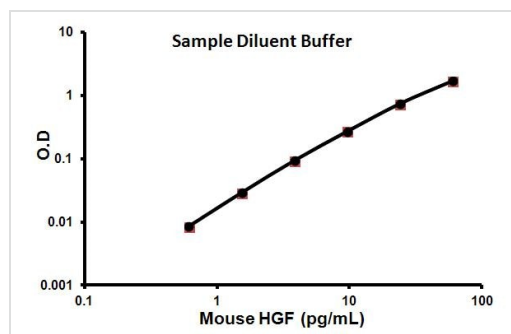
Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
2X Cell Lysis Buffer	1 x 5ml
500X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
5X Sample Diluent Buffer	1 x 10ml
Biotinylated anti-Mouse HGF	2 vials
HGF Microplate (12 x 8 wells)	1 unit
Recombinant Mouse HGF Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** Potent mitogen for mature parenchymal hepatocyte cells, seems to be a hepatotrophic factor, and acts as a growth factor for a broad spectrum of tissues and cell types. Activating ligand for the receptor tyrosine kinase MET by binding to it and promoting its dimerization.

**Involvement in disease** Deafness, autosomal recessive, 39

**Sequence similarities** Belongs to the peptidase S1 family. Plasminogen subfamily.  
Contains 4 kringle domains.  
Contains 1 PAN domain.  
Contains 1 peptidase S1 domain.

## Images



Representative Standard Curve using ab100687

Typical Standard Curve

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### **Our Abpromise to you: Quality guaranteed and expert technical support**

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors