

Human Met (c-Met) ELISA Kit ab277722

重组 SimpleStep ELISA

8 图像

概述

产品名称 人Met (c-Met) ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Serum	8			3.3%

批次间

样品	n	Mean	SD	CV%
Serum	3			4.4%

样品类型 Serum, Cell culture extracts, Cell Lysate, Cell culture media, Hep Plasma, EDTA Plasma

检测类型 Sandwich (quantitative)

灵敏度 81.5 pg/ml

范围 281.25 pg/ml - 18000 pg/ml

回收率 特定样本回收率

样品类型	平均%	范围
Serum	90	78% - 111%
Cell culture extracts	89	77% - 106%
Cell culture media	104	99% - 110%
Hep Plasma	104	91% - 113%
EDTA Plasma	103	99% - 108%
Cit plasma	94	84% - 113%

检测时间 1h 30m

实验步骤 One step assay

种属反应性 **与反应:** Human

产品概述 Human Met (c-Met) ELISA kit (ab277722) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Human Met (c-Met) protein in human cell and tissue extract samples. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Met (c-Met) with 81.6 pg/mL sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

说明 Met, also known as Hepatocyte Growth Factor Receptor (HGFR) or Proto-Oncogene c-Met (c-Met) is a receptor tyrosine kinase that transduces signals from the extracellular matrix into the cytoplasm by binding to hepatocyte growth factor ligand. Met regulates many physiological processes including proliferation, scattering, morphogenesis and survival. Ligand binding at the cell surface induces autophosphorylation of Met on its intracellular domain that provides docking sites for downstream signaling molecules. The recruitment of these downstream effectors by Met leads to the activation of several signaling cascades including the RAS-ERK, PI3 kinase-AKT, or PLC gamma-PKC. During embryonic development, Met signaling plays a role in gastrulation, development and migration of muscles and neuronal precursors, angiogenesis and kidney formation. In adults, Met participates in wound healing as well as organ regeneration and tissue remodeling. Met promotes also differentiation and proliferation of hematopoietic cells. Human and mouse Met are 90% identical on their amino acids sequences.

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.

平台 Pre-coated microplate (12 x 8 well strips)

性能

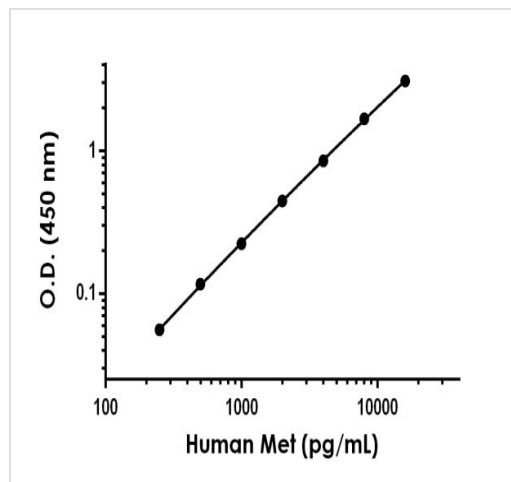
存放说明 Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Human Met (c-Met) Capture Antibody	1 x 600µl

组件	1 x 96 tests
10X Human Met (c-Met) Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 4BI	1 x 6ml
Human Met (c-Met) Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

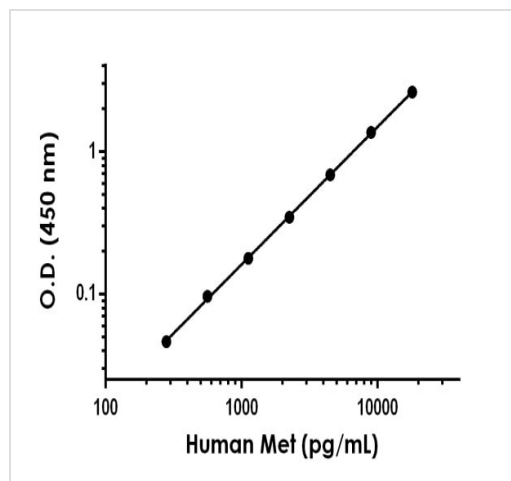
功能	Receptor for hepatocyte growth factor and scatter factor. Has a tyrosine-protein kinase activity. Functions in cell proliferation, scattering, morphogenesis and survival.
疾病相关	<p>Note=Activation of MET after rearrangement with the TPR gene produces an oncogenic protein.</p> <p>Note=Defects in MET may be associated with gastric cancer.</p> <p>Defects in MET are a cause of hepatocellular carcinoma (HCC) [MIM:114550].</p> <p>Defects in MET are a cause of renal cell carcinoma papillary (RCCP) [MIM:605074]. It is a subtype of renal cell carcinoma tending to show a tubulo-papillary architecture formed by numerous, irregular, finger-like projections of connective tissue. Renal cell carcinoma is a heterogeneous group of sporadic or hereditary carcinoma derived from cells of the proximal renal tubular epithelium. It is subclassified into common renal cell carcinoma (clear cell, non-papillary carcinoma), papillary renal cell carcinoma, chromophobe renal cell carcinoma, collecting duct carcinoma with medullary carcinoma of the kidney, and unclassified renal cell carcinoma.</p> <p>Note=A common allele in the promoter region of the MET shows genetic association with susceptibility to autism in some families. Functional assays indicate a decrease in MET promoter activity and altered binding of specific transcription factor complexes.</p> <p>Note=MET activating mutations may be involved in the development of a highly malignant, metastatic syndrome known as cancer of unknown primary origin (CUP) or primary occult malignancy. Systemic neoplastic spread is generally a late event in cancer progression. However, in some instances, distant dissemination arises at a very early stage, so that metastases reach clinical relevance before primary lesions. Sometimes, the primary lesions cannot be identified in spite of the progresses in the diagnosis of malignancies.</p>
序列相似性	<p>Belongs to the protein kinase superfamily. Tyr protein kinase family.</p> <p>Contains 3 IPT/TIG domains.</p> <p>Contains 1 protein kinase domain.</p> <p>Contains 1 Sema domain.</p>
结构域	The kinase domain is involved in SPSB1 binding.
翻译后修饰	Dephosphorylated by PTPRJ at Tyr-1349 and Tyr-1365.
细胞定位	Membrane.

图片



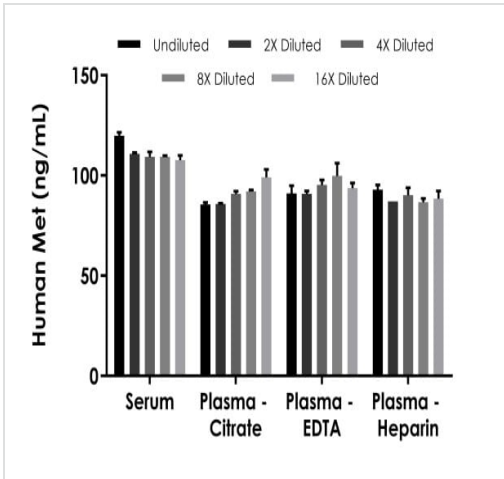
Example of human Met standard curve in Sample Diluent NS.

The Met standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



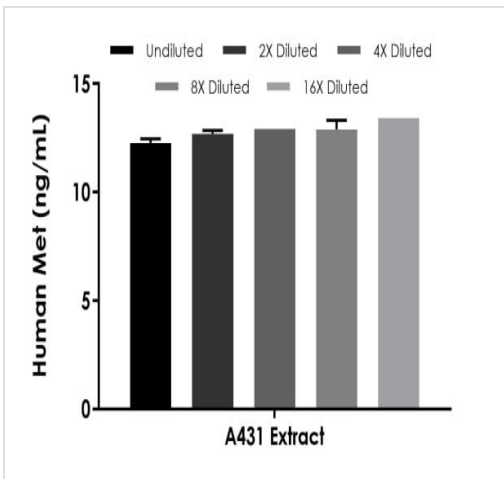
Extraction Buffer PTR.

The Met standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



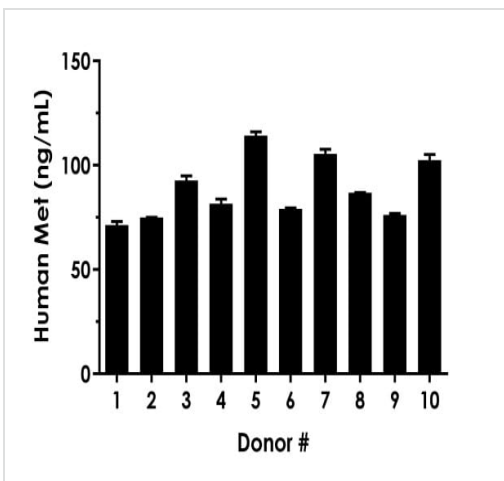
Interpolated concentrations of native Met in human serum and plasma samples.

The concentrations of Met were measured in duplicates, interpolated from the Met standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 10%, plasma (citrate) 10%, plasma (EDTA) 10% and plasma (heparin) 10%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Met concentration was determined to be 111.4 ng/mL in neat serum, 90.7 ng/mL in neat plasma (citrate), 94.1 ng/mL in neat plasma (EDTA) and 89.0 ng/mL in neat plasma (heparin).



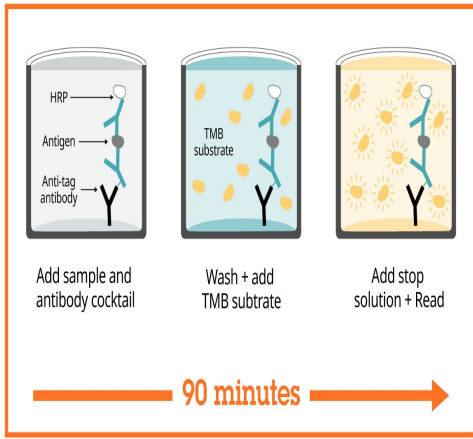
Interpolated concentrations of native Met in human A431 cell based on a 250 µg/mL extract load.

The concentrations of Met were measured in duplicate and interpolated from the Met standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Met concentration was determined to be 12.8 ng/mL in A431 extract.



Serum from ten individual healthy human female donors was measured in duplicate.

Interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Met concentration was determined to be 88.5 ng/mL with a range of 71.3 – 114.2 ng/mL.



Sandwich ELISA - Human Met (c-Met) ELISA Kit (ab277722)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Powered by recombinant antibodies

- Research with confidence**
Consistent and reproducible results
- Long-term and scalable supply**
Recombinant technology
- Success from the first experiment**
Confirmed specificity
- Ethical standards compliant**
Animal-free production

Recombinant Antibody Benefits

To learn more about the advantages of recombinant antibodies see [here](#).

Get more done with SimpleStep ELISA

- Easy to use**
Single-wash 90-minute protocol
- Flexible**
Matched antibody pairs available
- Precision antibodies**
High sensitivity, specificity and reproducibility
- Scalable**
Now in 10-pack and 384-well formats

Sandwich ELISA - Human Met (c-Met) ELISA Kit (ab277722)

To learn more about the advantages of SimpleStep ELISA[®] kits see [here](#).

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.cn/abpromise> or contact our technical team.

Terms and conditions

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