

Product datasheet

Mouse Tissue Factor Matched Antibody Pair Kit
ab218790

1 Image

Overview

Product name	Mouse Tissue Factor Matched Antibody Pair Kit
Detection method	Colorimetric
Assay type	ELISA set
Range	31.25 pg/ml - 2000 pg/ml
Species reactivity	Reacts with: Mouse
Product overview	Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant mouse Tissue Factor.

Both capture and detector antibodies are rabbit monoclonal antibodies.

Matched antibody pair kits and reagents deliver consistent, specific, and sensitive results.

Batch-to-batch consistency: only recombinant monoclonal antibodies are used in our matched antibody pairs.

Specificity: antibody pairs are screened in plasma and serum to ensure specificity in complex samples.

Sensitivity: benchmarked against commercially available antibody pairs to ensure equivalent or superior performance compared with the competition.

Additional buffers and plates are required for the assay. An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs ([ab210905](#)).

For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA® ([ab214091](#)) which uses the same antibody pair.

To receive an electronic copy of the Certificate of Analysis, please send an [email](#) with "CoA for matched antibody pair kit" in the subject line and the desired product number

and lot number in the body of the email.

Tested applications

Suitable for: ELISA

Platform

Reagents

Properties

Storage instructions

Store at -20°C. Please refer to protocols.

Components	10 x 96 tests	2 x 96 tests
Mouse Tissue Factor Capture Antibody	1 x 100µg	1 x 20µg
Mouse Tissue Factor Detector Antibody	1 x 25µg	1 x 5µg
Mouse Tissue Factor Lyophilized Protein	1 vial	1 vial

Function

Initiates blood coagulation by forming a complex with circulating factor VII or VIIa. The [TF:VIIa] complex activates factors IX or X by specific limited protolysis. TF plays a role in normal hemostasis by initiating the cell-surface assembly and propagation of the coagulation protease cascade.

Sequence similarities

Belongs to the tissue factor family.

Cellular localization

Membrane.

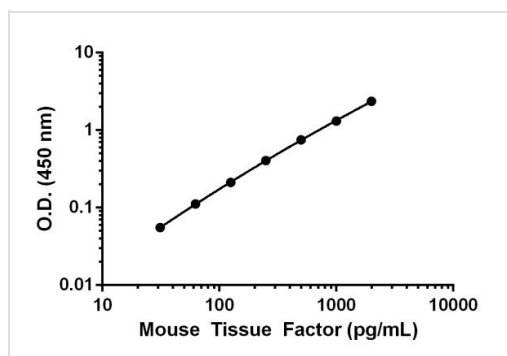
Applications

Our [Abpromise guarantee](#) covers the use of **ab218790** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.

Images



Mouse Tissue Factor standard curve

Standard calibration curve. Background subtracted values are graphed.

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