



ELISA PRODUCT INFORMATION & MANUAL

Human TFF3 ELISA Kit (Colorimetric) *NBP2-33318*

Enzyme-linked Immunosorbent Assay for quantitative detection. For research use only.

Not for diagnostic or therapeutic procedures.

NBP2-33318

Trefoil Factor 3, Intestinal (TFF3) ELISA Kit (Human)

Specificity

Human

Kit Type

Sandwich ELISA

Tests

4896

Sample Volume

100ul

Sensitivity

<23.8pg/ml

Detection Method

Colorimetric

Detection Range

62.5-4,000pg/ml

Sample Matrix

Serum, plasma, tissue homogenates, cell lysates, cell culture supernatants and other biological fluids

EU Commodity Code 38220000

UN DOT Shipping

UN2796 PGII

Shipping Temp

Blue Ice

Storage Temp

4°C/-20°C

The Trefoil Factor 3, Intestinal (TFF3) ELISA Kit (Human) is a sandwich enzyme immunoassay for the in vitro quantitative measurement of TFF3 in human serum, plasma, tissue homogenates, cell lysates, cell culture supernatants and other biological fluids.

Detection Range

62.5-4,000pg/ml

Sensitivity

<23.8pg/ml

Precision

Intra-Assay: CV<10% Inter-Assay: CV<12%

Test Principle

The microtiter plate provided in this kit has been pre-coated with an antibody specific to TFF3. Standards or samples are then added to the appropriate microtiter plate wells with a biotin-conjugated antibody specific to TFF3. Next, Avidin conjugated to Horseradish Peroxidase (HRP) is added to each microplate well and incubated. After TMB substrate solution is added, only those wells that contain TFF3, biotin-conjugated antibody and enzyme-conjugated Avidin will exhibit a change in color.

The enzyme-substrate reaction is terminated by the addition of sulfuric acid solution and the color change is measured spectrophotometrically at a wavelength of 450nm ± 10nm. The concentration of TFF3 in the sample is then determined by comparing the O.D. of the sample to the standard curve.

Kit Components

*Microtiter Plate, 1x96 wells. Pre-coated; ready to use.

*Standard, 2x1vial

*Standard Diluent, 1x20ml

*Detection Reagent A, 1x120ul

*Detection Reagent B, 1x120ul

*Assay Diluent A, 1x12ml

*Assay Diluent B, 1x12ml

*TMB Substrate, 1x9ml

*Stop Solution, 1x6ml

*Wash Buffer, 30X, 1x20ml

Storage and Stability

Store Microtiter Plate, Standard, Detection Reagent A, and Detection Reagent B at -20°C. Store all the other components at 4°C. Unused kit is stable for 6 months. Once kit components are opened, it is highly recommended to use remaining reagents within 1 month provided this is within the expiration date of the kit. For maximum recovery of product, centrifuge the original vials after thawing and prior to removing the cap.

Assay Procedure Summary

1. Prepare all reagents, samples and standards. 2. Add 100ul standard or sample to each well. Incubate 1 hour at 37°C. 3. Aspirate and add 100ul Detection Reagent A. Incubate 1 hour at 37°C. 4. Aspirate and wash 3 times. 5. Add 100ul Detection Reagent B. Incubate 30 minutes at 37°C. 6. Aspirate and wash 5 times. 7. Add 90ul TMB Substrate. Incubate 10-20 minutes at 37°C. 8. Add 50ul Stop Solution. Read at 450nm immediately.

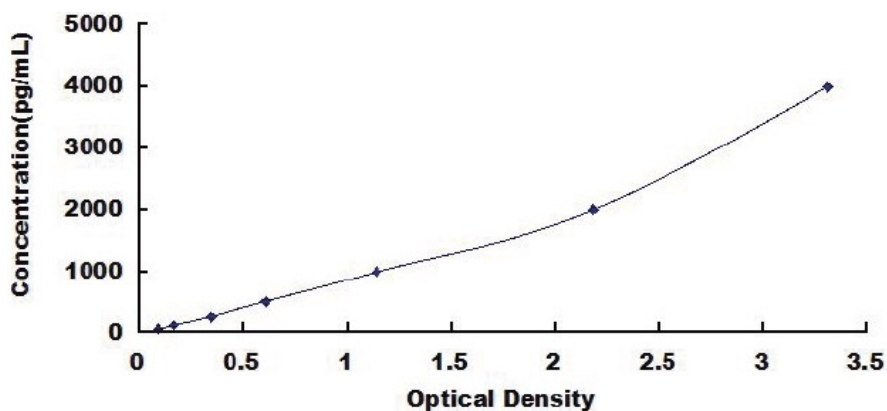


Fig. 1 Typical standard curve for Trefoil Factor 3, Intestinal (TFF3) ELISA Kit (Human) (shown for reference only; user must generate a new standard curve with each assay).