

# Human MIP3 $\alpha$ ELISA kit (DDXK-E-MIP3 $\alpha$ )

#### Intended use

The human MIP3 $\alpha$  ELISA kit is for quantitative determination of MIP3 $\alpha$  concentrations in human cell culture supernatants and biological fluids (plasma, serum). This kit is for research purpose only.

## **Background**

Macrophage inflammatory protein (MIP3α)/ CCL20 is a CC-type chemokine mapped to chromosome 2 in humans. Langerhans cells (LCs) represent a unique population of DCs colonizing epithelium. MIP3 $\alpha$  plays a central role in LC precursor recruitment into the epithelium during inflammation. (a) Among DC populations, MIP3 $\alpha$  is the most potent chemokine inducing the selective migration of in vitro-generated CD34<sup>+</sup> hematopoietic progenitor cell-derived LC precursors and skin LCs, in accordance with the restricted MIP3 $\alpha$  receptor (CC chemokine receptor 6) expression by these cells. (b) MIP3 $\alpha$  is mainly produced by epithelial cells, and the migration of LC precursors induced by the supernatant of activated skin keratinocytes is blocked by an antibody against MIP3α. (c) In vivo, MIP3 $\alpha$  is selectively produced at sites of inflammation as illustrated in tonsils and lesional psoriatic skin where MIP3 $\alpha$  up regulation is associated with an increase in LC turnover. (d) The secretion of MIP3 $\alpha$  is strongly up regulated by cells of epithelial origin after inflammatory stimuli (interleukin 1 $\beta$ plus TNFα) or T cell signals (Homey B. et al, 2000, J.Immunol. 164:6621-32). Monoclonal antibodies anti-MIP3α were obtained after mice immunization with MIP3α-transfected eukaryotic cells. These clones were used to set up a human MIP3 $\alpha$  specific Elisa, which also detects macaque MIP3 $\alpha$ . No cross-reactivity with MIP1α, MIP1β, MIP1β, MIP3β, eotaxin, 6Ckine, Rantes, MCP-1, TARC, MDC, TECK, SDF-1 and IP-10 was detected (Dieu-Nosjean et al, J;Exp.med, 2000 192: 705-18).

#### **Kit contents**

*Capture Antibody*: 0.5mg/mL of mouse anti-MIP3 $\alpha$  monoclonal antibody (#DDX0430, clone 319F6).

**Detection Antibody**: 0.5mg/mL of HRPO-conjugated anti-MIP3 $\alpha$  mouse monoclonal antibody (#DDX0420, clone 206D9).

**Standard**: Each vial contains  $1\mu g/mL$  of recombinant human MIP3 $\alpha$  produced, purified and concentrated from eukaryotic cells. A 7-point standard curve using 2-fold serial dilutions in sample dilution buffer, and a high standard of 20 ng/mL is recommended.

#### Storage

All the reagents should be aliquoted before storage. Minimize repeated freeze and thaw. Refer to expiration date on all reagents prior to use. Do not mix reagents from different kits unless they have the same lot number. Capture and detection antibodies storage: -20°C

## Materials and reagents required but not provided

96well-plate Nunc Maxi Sorp 50mM Carbonate buffer pH9.6 PBS-1% BSA-0.05% Tween20 TMB super sensitive HRP (TMBS100-0500, TEBU-BIO) Multichannel pipettes and pipette tips A standard microplate reader (620nm)



### Sensitivity

The minimum detectable dose of human MIP3 $\alpha$  was determined to be approximately 0.2ng/ml.

#### Principle of the assay

The human MIP3 $\alpha$  ELISA kit is for the quantitative determination of human MIP3 $\alpha$  in human cell culture supernatants, plasma, serum, and various biological fluids. This ELISA kit contains the specific components required for the development of human MIP3 $\alpha$  sandwich ELISAs. Each kit contains sufficient materials to run ELISAs on 3X96-well plates.

The DDX MIP3 $\alpha$  ELISA kit is a solid phase sandwich ELISA (Enzyme-Linked Immunosorbent Assay). The capture monoclonal antibody specific for MIP3 $\alpha$  is coated on a 96-well plate. Standards and samples are added to the wells, and any MIP3 $\alpha$  present binds to the immobilized antibody. The wells are washed and a HRPO-conjugated anti-MIP3 $\alpha$  monoclonal antibody is then added, producing an antibody-antigen-antibody "sandwich". To produce color in proportion to the amount of MIP3 $\alpha$  present in the sample, TMB substrate solution is loaded and absorbance is measured at 620 nm.

#### **Plate Preparation**

## Pre-warm all the reagents to room temperature prior to setting up the assay

- 1. Dilute the capture antibody to  $2.5\mu g/ml$  in 50mM carbonate buffer pH9.6. Immediately coat a 96-well microplate with  $120\mu L$  per well of the diluted capture antibody. Seal the plate and incubate overnight at room temperature (RT).
- 2. Remove coating solution, and wash with at least 200µl of PBS-0.05% Tween20, repeating the process two times for a total of three washes. Complete removal of liquid at each step is essential for good performance. After the last wash, remove any remaining wash buffer by inverting the plate and blotting it against clean paper towels.

The plates are now ready for sample addition.

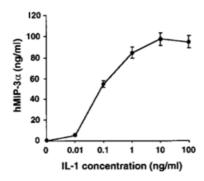
## **Assay Procedure**

- 1. Add  $100\mu$ L/well of sample or standards, diluted in PBS-1% BSA-0.05% Tween20. Seal the plate and incubate 1h30 at 37°C.
- 2. Repeat the aspiration/wash as in step 2 of plate preparation.
- 3. Add  $100\mu$ L/well of the HRPO-conjugated detection antibody, diluted in PBS-1% BSA-0.05% Tween20 to  $3\mu$ g/ml. Seal the plate and incubate 1h30 at 37°C.
- 4. Repeat the aspiration/wash as in step 2 of plate preparation.
- 5. Add 100μL/well of TMB. Incubate for 5-10 minutes at RT.
- 6. Determine the optical density of each well, using a microplate reader set to 620 nm.



#### Standard curve

Each laboratory should establish its own standard curve. Here is an example of MIP3 $\alpha$  dosage using DDX-K-E-MIP3 $\alpha$ .



Elisa analysis of MIP-3lpha secretion by CHA cell line cultured in the presence of increasing doses of IL1-eta

Dieu-Nosjean et al, 2000

## **Troubleshootings**

To obtain good and reproducible results, usage of sterile reagents and clean materials is strongly recommended. All basic reagents such as washing and dilution buffers, water, must be devoid of contamination.

To ensure pH stability, incubation at  $37^{\circ}\text{C}$  should be performed in a humidified atmosphere of 5%  $\text{CO}_2$ 

Problems	Possible Sources	Solutions
Poor detection value	The concentration of antigen in samples was too low	Enriching samples to increase the concentration
	Samples were ineffective	Check if the samples are stored at cold
		environment. Detect samples in timely manner
No signal	Incorrect or no Detection Antibody was added	Add appropriate Detection Antibody and continue
	Substrate solution was not added	Add substrate solution and continue
	Incorrect storage condition	Check if the kit is stored at recommended
		condition and used before expiration date
Poor Standard Curve	Inappropriate storage	Aliquot standard and store at -70°C
	Imprecise / inaccurate pipetting	Check / calibrate pipettes
	Incubations done at inappropriate temperature or timing	Follow the general ELISA protocol
	Background wells were contaminated	Avoid cross contamination by using the sealer
High Background	Insufficient washes	Use multichannel pipettes without touching the
		reagents on the plate
		Increase cycles of washes and soaking time
	TMB Substrate Solution was contaminated	TMB Substrate Solution should be clear and
		colorless prior to addition to wells
	Materials were contaminated	Use clean plates, tubes and pipettes tips
Non-specificity	Samples were contaminated	Avoid cross contamination of samples
	The concentration of samples was too high	Try higher dilution rate of samples