### **5**-

# Hydroxymethylcytosine DNA ELISA kit

## Highly sensitive, rapid ELISA kit to quantify 5-hydroxymethylcytosine in DNA with results

That-Hydroxymethyleytopine PNATASA sipile a complete, golorimetric, immunometric immunoassay kit for the quantitative determination of 5-analysis, hydroxymethylcytosine in DNA with results in just 3 hours.

The 5-methylcytosine (5-mC) modification of DNA in epigenetic regulation has been well studied over the last several decades. However, the role of the so-called "sixth base", 5-hydroxymethylcytosine (5-hmC), has yet to be defined. Recent studies have associated 5-hmC patterns to transcriptional regulation of genes, which may contribute to normal and disease states of organisms.

Citations: 2

View Online »

**Ordering Information** 

Order Online »

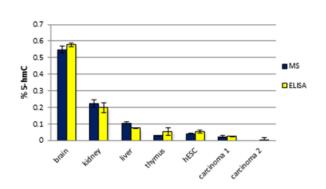
ADI-900-225-0001

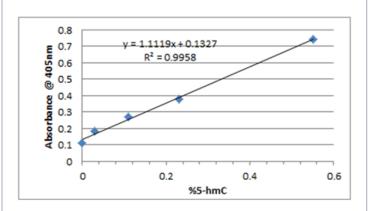
96 wells

Manuals, SDS & CofA

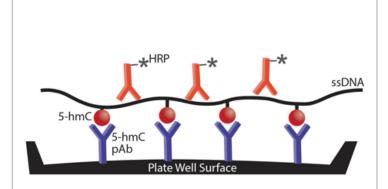
View Online »

- Rapid fully quantitative results in
  3 hours
- Convenient user-friendly alternative to mass spectrometry
- Simple efficient and wellestablished protocol
- High-throughput 96 well format ideal for global detection, tissuespecific quantitation, highthroughput compound screening, and more





DNA Mass Spectroscopy versus ELISA analysis: The 5-hmC DNA ELISA Kit (prod. no. ADI-900-225) quantifies 5-hmC in numerous DNA samples with close correlation to LC-MS/MS-MRM analysis. DNA samples include: brain, kidney, liver, thymus, human embryonic stem cells and two carcinoma samples.



#### **Handling & Storage**

**Use/Stability** Store all components at 4°C, except antibody and control set should be stored at -20°C,

while the DNA HRP antibody should be stored at -80°C.

Shipping Dry Ice

#### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name 5-hmC

**Application** Colorimetric detection, ELISA

**Application Notes**For the quantitative determination of species independent

5-hydroxymethylcytosine DNA in a variety of DNA

samples.

Assay Time ~3 hours

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

Contents Microtiter Plate, Coating Buffer, ELISA Buffer, Primary

Antibody, Antibody-conjugate, HRP Developer, Controls

Sensitivity <0.02% 5-hydroxymethylcytosine DNA per 100ng input

DNA (range 0.03-0.55% (100 ng/µl))

Species Reactivity Species independent

Wavelength 450 nm

Last modified: May 29, 2024

