

# 5-

## Hydroxymethylcytosine DNA ELISA kit

### Highly sensitive, rapid ELISA kit to quantify 5-hydroxymethylcytosine in DNA with results that closely correlate to mass spectrometry analysis.

The 5-Hydroxymethylcytosine DNA ELISA kit is a complete, colorimetric, immunometric immunoassay kit for the quantitative determination of 5-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 »](#)

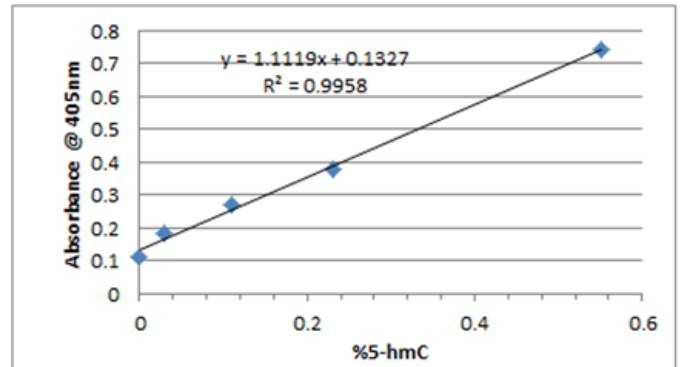
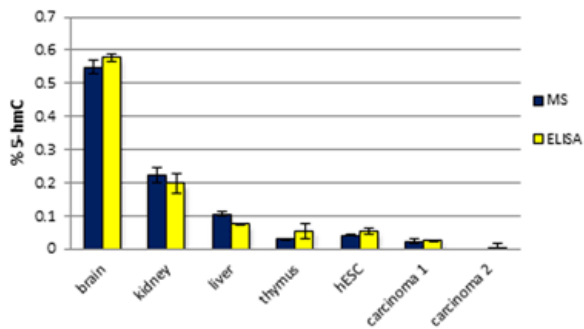
ADI-900-225-0001

96 wells

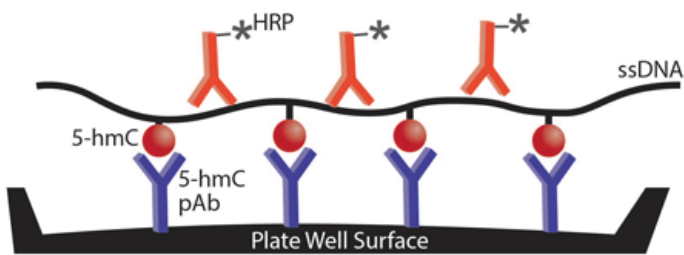
### Manuals, SDS & CofA

[View Online »](#)

- Rapid - fully quantitative results in < 3 hours
- Convenient – user-friendly alternative to mass spectrometry
- Simple – efficient and well-established protocol
- High-throughput – 96 well format ideal for global detection, tissue-specific quantitation, high-throughput 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

**Species Reactivity** Species independent

**Wavelength** 450 nm

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)