

Global 5-hmC DNA ELISA Kit

Catalog No: 55025

Format: 1 x 96 rxns

Quality Control: The Global 5-hmC Quantification Kit is tested for sensitivity in detecting changes in global 5-hydroxymethylcytosine (5-hmC) levels in genomic DNA. 96-stripwell plates are coated with a DNA binding agent and *Mse*I digested genomic DNA or serial dilutions of the included 5-hmC DNA Standard synthetic oligo are added to each well. A primary antibody specific for 5-hmC and a secondary antibody conjugated to HRP are used for detection. The colorimetric readout is quantified by spectrophotometry using a microplate reader at 450 nm.

The OD 450 nm results are graphed using Excel Linear Regression. The standard curve for Lot 1272001 with a 6 minute developing time is shown in Figure 1. The ng 5-hmC and % 5-hmC are calculated for the DNA samples from the standard curve using the trendline equation.

Storage and Guarantee: Components can be stored at -20°C prior to first use, then we recommend storing the components at the temperatures listed in the manual. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

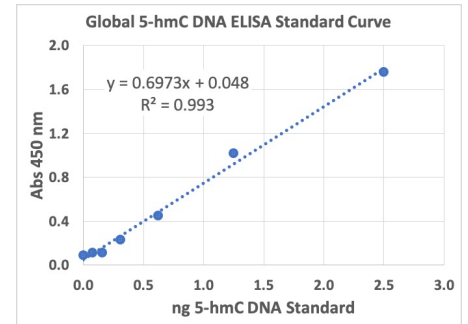


Figure 1: Global 5-hmC DNA ELISA Standard Curve