

**Amplite® Colorimetric Beta-Hydroxybutyrate (Ketone Body) Assay Kit**

Catalog number: 13830

Unit size: 200 Tests

Component	Storage	Amount (Cat No. 13830)
Component A: Enzyme Mix	Freeze (< -15 °C), Minimize light exposure	2 bottles (lyophilized powder)
Component B: Assay Buffer	Freeze (< -15 °C)	1 bottle (10 mL)
Component C: NAD	Freeze (< -15 °C), Minimize light exposure	1 vial
Component D: $\beta$ -Hydroxybutyrate ( $\beta$ -HB) Standard	Freeze (< -15 °C), Minimize light exposure	1 vial (10 $\mu$ L)

**OVERVIEW**

Ketone bodies are produced by the liver and used peripherally as an energy source when blood glucose levels drop. The two main ketone bodies are beta-hydroxybutyrate (beta-HB) and acetoacetate (AcAc), while acetone is the third abundant ketone body. Normally these two predominant ketone bodies are present in small amounts in the blood during fasting (low food intake) and prolonged exercise. In patients who have diabetes, alcohol or salicylate poisoning, hormone deficiency, childhood hypoglycemia and other acute disease states, large quantities of ketone bodies are found in the blood. The over-production and accumulation of ketone bodies in the blood (ketosis) can lead to pathological metabolic acidosis (ketoacidosis). In extreme cases, ketoacidosis can be fatal. Blood ketone testing methods that quantify beta-HB, the predominant ketone body in the blood (approximately 75%) have been used for diagnosing and monitoring treatment of ketoacidosis. Amplite® Colorimetric Beta-Hydroxybutyrate Assay Kit offers a sensitive colorimetric assay for measuring beta-HB levels in biological samples. This assay is based on an enzyme coupled reaction of beta-HB, in which the NADH product can be specifically monitored by a NADH sensor. The signal can be measured by an absorbance microplate reader with the OD ratio at the wavelength of 570 nm to 610 nm. With this Amplite® Colorimetric Beta-Hydroxybutyrate Assay Kit, we were able to detect as low as 4  $\mu$ M beta-HB in a 100  $\mu$ L reaction volume.

**AT A GLANCE**
**Protocol Summary**

1. Prepare  $\beta$ -HB working solution (50  $\mu$ L)
2. Add  $\beta$ -HB standards or test samples (50  $\mu$ L)
3. Incubate at room temperature for 10 - 30 min
4. Monitor Absorbance increase at OD ratio of 570/610 nm

**Important Note**

Thaw one vial of each kit component at room temperature before starting the experiment.

**KEY PARAMETERS**
**Absorbance microplate reader**

Absorbance	570/610 nm
Recommended plate	Clear bottom

**PREPARATION OF STOCK SOLUTIONS**

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles

**NAD stock solution (100X)**

Add 100  $\mu$ L of H<sub>2</sub>O into the vial of NAD (Component C) to make 100X NAD stock solution.

 **$\beta$ -HB standard solution (100 mM)**

Add 1 mL of H<sub>2</sub>O or 1X PBS buffer into the vial of  $\beta$ -HB standard (Component D) to make 100 mM  $\beta$ -HB standard solution.

**PREPARATION OF STANDARD SOLUTIONS**

For convenience, use the Serial Dilution Planner:  
<https://www.aatbio.com/tools/serial-dilution/13830>

 **$\beta$ -HB standard**

Add 10  $\mu$ L of 100 mM  $\beta$ -HB standard stock solution into 990  $\mu$ L 1X PBS buffer to generate 1000  $\mu$ M  $\beta$ -HB standard solution (HB7). Take the 1000  $\mu$ M  $\beta$ -HB standard solution (HB7) and perform 1:3 serial dilutions in PBS to get serially diluted  $\beta$ -HB standards (HB6 - HB1). Note: Diluted  $\beta$ -HB standard solution is unstable, and should be used within 4 hours.

**PREPARATION OF WORKING SOLUTION**

Add 50  $\mu$ L of 100X NAD stock solution into the bottle of Component A+B, and mix well to make  $\beta$ -HB working solution (Component A+B+C).

Add 5 mL of Assay Buffer (Component B) into one bottle of Enzyme Mix (Component A).

**Note:** This  $\beta$ -HB working solution is not stable, use it promptly and avoid direct exposure to light.

**SAMPLE EXPERIMENTAL PROTOCOL**

**Table 1.** Layout of  $\beta$ -HB standards and test samples in a clear bottom 96-well microplate. HB =  $\beta$ -HB standard (HB1 - HB7, 1 to 1000  $\mu$ M); BL = blank control; TS = test sample.

BL	BL	TS	TS
HB1	HB1	...	...
HB2	HB2	...	...
HB3	HB3		
HB4	HB4		
HB5	HB5		
HB6	HB6		
HB7	HB7		

**Table 2.** Reagent composition for each well.

Well	Volume	Reagent
HB1-HB7	50 $\mu$ L	Serial Dilution (1 to 1000 $\mu$ M)
BL	50 $\mu$ L	1X PBS Buffer
TS	50 $\mu$ L	Test Sample

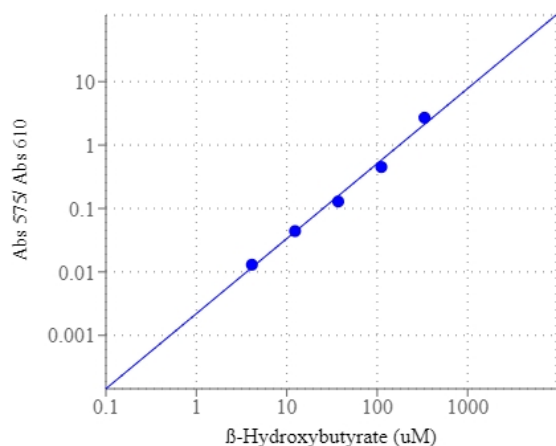
1. Prepare  $\beta$ -HB standards (HB), blank control (BL) and test samples (TS) according to the layout provided in Tables 1 and 2. For a 384-well plate, use 25  $\mu$ L of reagent per well instead of 50  $\mu$ L.
2. Add 50  $\mu$ L of  $\beta$ -HB working solution to each well of  $\beta$ -HB standard, blank control, and test samples to make the total assay volume of 100  $\mu$ L/well. For a 384-well plate, add 25  $\mu$ L of  $\beta$ -HB working solution into each well instead, for a total volume of 50  $\mu$ L/well.
3. Incubate the reaction at room temperature for 10 - 30 minutes, protected from light.
4. Monitor the absorbance increase with an absorbance plate reader at OD ratio of 570/610 nm.

#### EXAMPLE DATA ANALYSIS AND FIGURES

The reading (Abs 575/ Abs 610) obtained from the blank standard well is used as a negative control. Subtract this value from the other standards' readings to obtain the base-line corrected values. Then, plot the standards' readings to obtain a standard curve and equation. This equation can be used to calculate  $\beta$ -Hydroxybutyrate samples. We recommend using the Online Linear Regression Calculator which can be found at:

<https://www.aatbio.com/tools/linear-logarithmic-semi-log-regression-online-calculator>

Chart



**Figure 1.**  $\beta$ -Hydroxybutyrate ( $\beta$ -HB) dose response was measured with Amplite™ Colorimetric  $\beta$ -Hydroxybutyrate Assay Kit on a black wall/clear bottom 96-well plate using a SpectraMax microplate reader (Molecular Devices).

#### DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.