

# **Amplite® Rapid Colorimetric Biotin Quantitation Kit \*Optimized to Use with Nanodrop\***

Catalog number: 5537  
Unit size: 100 tests

Component	Storage	Amount (Cat No. 5537)
Component A: Avidin	Freeze (< -15 °C), Minimize light exposure	1 Vial
Component B: HABA Assay Buffer	Freeze (< -15 °C), Minimize light exposure	2 Vials (1 mL/Vial)
Component C: d-Biotin	Freeze (< -15 °C), Minimize light exposure	1 Vial (200 µL, 100 µM)

## **OVERVIEW**

Amplite™ Colorimetric Biotin Quantitation Kit is designed to quantify the number of biotins attached to a protein or other macromolecule with a Nanodrop Spectrometer. The kit uses HABA (4'-hydroxyazobenzene-2-carboxylic acid) reagent that demonstrates a dramatic spectral change when bound to avidin. Biotin easily displaces HABA from the HABA/avidin complex, resulting in a decrease of absorption at 500 nm. By measuring the absorbance change before and after the addition of the biotin-containing sample at 500 nm, the amount of biotin in a sample can be quantified spectrophotometrically. The kit is best used to determine biotin concentration in the range from 2 to 16 µM. It is optimized to use with Nanodrop spectrometer. Biotin is a relatively small molecule that is routinely conjugated to antibodies and proteins with minimal interference of their biological activity. The avidin/streptavidin-biotin interaction is the strongest known binding pair between a protein and its ligand. The biotin-avidin interaction has been extensively explored for a variety of biological applications.

## **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*

### **100X Avidin Stock Solution**

1. To prepare a 100X Avidin stock solution, add 40 µL of ddH<sub>2</sub>O into the vial of Avidin (Component A), and mix thoroughly.

**Note:** Any unused 100X Avidin stock solution should be divided into single-use aliquots and stored at -20°C.

## **PREPARATION OF WORKING SOLUTION**

### **HABA/Avidin Working Solution**

1. To prepare a HABA/avidin working solution, add 2 µL of the 100X Avidin stock solution into 200 µL of the HABA assay buffer (Component B), and mix thoroughly.

**Note:** Any unused HABA/avidin working solution can be store for up to one week at 4°C.

## **SAMPLE EXPERIMENTAL PROTOCOL**

### **Quantitate Amount Biotin with a Nanodrop**

1. Prepare the following controls and samples as outlined in the

table below:

Positive Control	Negative Control	Test Sample*
d-Biotin (Component C): 2 µL	ddH <sub>2</sub> O: 2 µL	2 µL
HABA/Avidin working solution: 18 µL	HABA/Avidin working solution: 18 µL	HABA/Avidin working solution: 18 µL

**Note:** To ensure the concentration of biotin falls within the assay's linear range (2-16 µM of biotin final concentration), it is necessary to test the biotin-containing samples at several dilutions.

**Note:** To ensure assay accuracy, avoid using buffers containing potassium, as it causes precipitation.

**Note:** Free biotin must be separated from the biotinylated protein by gel filtration or dialysis.

2. Incubate the reaction mixture at room temperature for 5 minutes, protected from light. Avoid creating bubbles during pipetting.
3. Monitor absorbance at 500 nm using a Nanodrop Spectrometer, with a sample measurement of 2~2.5 µL. Utilize the protein label function for readings.

## **DATA ANALYSIS - CALCULATIONS**

### **1. Calculate the change of absorbance at 500 nm:**

$$\Delta A_{500} = A_{500} \text{ of negative control} - A_{500} \text{ of Biotin sample or positive control}$$

### **2. Calculate the biotin concentration (M):**

$$M = [\Delta A_{500} \div \epsilon_{\text{HABA/Biotin}}] \times \text{DF}$$

Where:

- $\epsilon_{\text{HABA/Biotin}} = 34,500 \text{ M}^{-1}\text{cm}^{-1}$
- DF = Dilution Factor

### **3. Calculate protein concentration (P)**

$$P = \text{protein concentration (mg/mL)} \div \text{molecular weight of protein}$$

### **4. Calculate molar ratio of biotin to protein (MR)**

$$\text{MR} = M/P$$

Where:

- M = Biotin Concentration
- P = Protein Concentration

#### 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.