

HABA Biotin Quantification Kit *Colorimetric*

<i>Revision#1.3</i>	<i>Last Updated: September 2021</i>
Catalog #	AS-72096-120
Kit Size	25 Cuvette Assays or 120 Microplate Assays

- **Optimized Performance:** Optimal conditions for the quantification of biotin.
- **Enhanced Value:** Less expensive than the sum of individual components.
- **High Speed:** Minimal hands-on time.
- **Assured Reliability:** Detailed protocol and references are provided.

Kit Components, Storage and Handling

Component	Description	Quantity
Component A	Avidin	1 bottle
Component B	HABA assay buffer	25 mL
Component C	d-biotin	100 µM, 200 µL

Other Materials Required (but not provided)

- 96-well microplate or cuvettes: Clear microplate or cuvettes.
- Absorbance microplate reader or spectrophotometer: Capable of detecting absorbance at 500 nm.

Storage and Handling

- Store Component A at -20 °C.
- Store all other components at 4°C

Introduction

Biotin and avidin (or streptavidin) bind non-covalently with a higher binding affinity than most antigen-antibody interactions. This very tight binding makes labeling proteins with biotin a useful tool for applications such as affinity chromatography and immunoanalytical methods. Reaction conditions for biotinylation are chosen such that the target molecule (e.g. an antibody) is labeled with enough biotin residues to purify or detect the molecule, but not so much that the biotin interferes with the function of the molecule.

The HABA Biotin Quantitation Kit provides a convenient method to estimate the molar ratio of biotin to protein on biotinylated conjugates. It also can be used to quantitate the biotin concentration in a solution. The assay utilizes the observation that HABA (4'-hydroxyazobenzene-2-carboxylic acid) shows dramatic spectral changes when it binds to avidin. Free HABA has an absorption peak at 348 nm, while the HABA/avidin complex has strong absorption at 500 nm. Since the affinity between HABA and avidin is relatively weak ($K_d=5.8 \times 10^{-6}$ M) compared to the affinity between biotin and avidin ($K_d=1 \times 10^{-15}$ M), biotin can easily replace HABA from the HABA/avidin complex, resulting in a decrease of absorption at 500 nm.

Avidin and HABA at an optimal ratio are included in the assay kit. The kit has a linear range 2 to 16 μ M of biotin (final concentration). Assay can be performed in cuvette or microplate format.

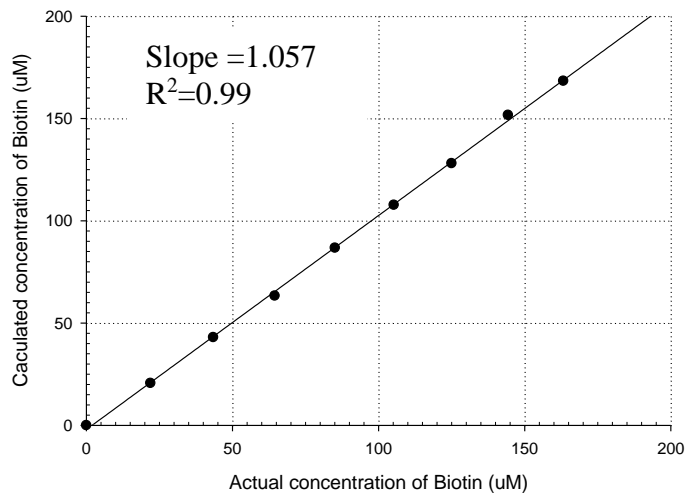


Figure 1. The calculated concentration of biotin is consistent with the actual concentration of biotin.

Protocol

Protocol A Quantitating Biotin in a Microplate Format

Note 1: It is necessary to test the biotin-containing sample at several dilutions to ensure that the concentration of biotin is within the assay linear range, 2-16 μ M of biotin (final concentration). You can roughly estimate the biotin concentration using the protein concentration of your conjugate solution. For example, biotinylated IgG at concentration 1 mg/mL contains 6.6 μ M biotin if the labeling ratio is 1 biotin per 1 IgG molecule. If the conjugate is too diluted, you will need to concentrate the solution in order to accurately measure biotin concentration. Protein solution can be concentrated by using a speed vacuum or a centrifugal filter (Millipore, Cat# MRCPRT010).

Note 2: Avoid buffers containing potassium, as it will cause precipitation in the assay.

Note 3: Free biotin must be separated from the biotinylated protein by dialysis or gel filtration.

1. Prepare HABA/Avidin assay mixture.

- Add 25 mL of HABA assay buffer (Component B) into one bottle of avidin (Component A). Mix the reagent completely.

Note: The unused portion of HABA/Avidin assay mixture may be stored at 4°C up to one week.

2. Biotin Assay:

- Add 180 μ L of HABA/Avidin assay mixture per well in a 96-well plate. Besides the wells for test samples, prepare extra wells for negative and positive controls.
- Add 20 μ L of biotin-containing sample into each well.

- Add 20 μL of deionized water or the same buffer used to dissolve biotin-containing sample into the negative control wells.
- Add 20 μL of d-biotin (Component C) into the positive control wells.
- Mix the reagents well by shaking on a plate shaker at 100-200 rpm for 5 min. Avoid creating bubbles during pipetting.
- Read absorbance at 500 nm.

3. **Data analysis:** Refer to the Data Analysis Section.

Protocol B Quantitating Biotin in a Cuvette Format

Note 1: It is necessary to test the biotin-containing sample at several dilutions to ensure that the concentration of biotin is within the assay linear range, 2-16 μM of biotin (final concentration). You can roughly estimate the biotin concentration through the protein concentration of your conjugate solution. For example, 1 mg/mL biotinylated IgG contains 6.6 μM biotin if the labeling ratio is 1 biotin per IgG molecule. If the conjugate is too diluted, you will need to concentrate the solution in order to accurately measure biotin concentration. Protein solution can be concentrated by using a speed vacuum or a centrifugal filter (Millipore, Cat# MRCPRT010).

Note 2: Avoid buffers containing potassium, as it will cause precipitation in the assay.

Note 3: Free biotin must be separated from the biotinylated protein by dialysis or gel filtration.

1. Prepare HABA/Avidin assay mixture.

- Add 25 mL of HABA assay buffer (Component B) into one bottle of avidin (Component A). Mix the reagent completely.

Note: The unused portion of HABA/Avidin assay mixture may be stored at 4°C up to one week.

2. Biotin Assay:

- Add 900 μL of HABA/Avidin assay mixture into a cuvette. Read absorbance at 500 nm. Record this reading as $A_{500\text{nm}}$ of HABA/Avidin negative control.
- Positive control: Add 900 μL of HABA/Avidin assay mixture into a cuvette. Add 100 μL of d-biotin (Component C) into the cuvette. Mix the reagents thoroughly and avoid creating bubbles. Read absorbance at 500 nm. Record the reading as $A_{500\text{nm}}$ of HABA/Avidin/Biotin sample for the positive control.
- Add 100 μL of biotin-containing sample into the cuvette. Mix the reagents well and avoid creating bubbles. Read absorbance at 500 nm. Record the reading as $A_{500\text{nm}}$ of HABA/Avidin/Biotin sample.

3. **Data analysis:** Refer to the Data Analysis Section.

Data Analysis

Calculate data from a microplate format

1. $\Delta A_{500\text{nm}} = A_{500\text{nm}}$ of negative control - $A_{500\text{nm}}$ of Biotin sample or positive control
2. Biotin concentration (M) = $[\Delta A_{500\text{nm}} / (34,500 \times 0.5)] \times \text{dilution factor}$
Note: $\epsilon_{\text{HABA/Avidin}} = 34,500 \text{ M}^{-1}\text{cm}^{-1}$, Light path = 0.5 cm
3. Protein concentration (M) = protein concentration (mg/mL) / molecular weight of protein
4. Molar ratio of biotin to protein = Biotin concentration (M) / Protein Concentration (M)

Calculate data from a cuvette format

1. $A_{500\text{ nm}} = (0.9 \times A_{500\text{ nm of HABA/Avidin}}) - A_{500\text{ nm of HABA/Avidin/Biotin sample}}$
2. Biotin concentration (M) = $(\Delta A_{500\text{ nm}} / 34,500) \times \text{dilution factor}$
Note: $\epsilon_{\text{HABA/Avidin}} = 34,500 \text{ M}^{-1}\text{cm}^{-1}$, Light path = 1 cm
3. Protein concentration (M) = protein concentration (mg/mL) / molecular weight of protein
4. Molar ratio of biotin to protein = Biotin concentration (M) / Protein Concentration (M)

Positive control

1. Calculate the biotin concentration of the positive control according to the above formulas and then compare it to its actual concentration of 100 μM . The calculated concentration should be very close to the actual concentration if all the procedures and calculations have been correctly performed.