

Product Information

Ketone Body Assay Kit

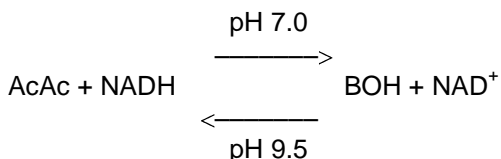
Catalog Number **MAK134**Storage Temperature -20°C

TECHNICAL BULLETIN

Product Description

Ketone bodies, 3-hydroxybutyric acid (BOH) and acetoacetic acid (AcAc), are produced in the liver primarily from oxidation of fatty acids, and are normally present at low concentrations in urine and blood. Increased ketone concentrations in the blood may lead to metabolic acidosis, which has been associated with diabetes, childhood hypoglycemia, growth hormone deficiency, alcohol or salicylate intoxication, and inborn errors of metabolism.

In this kit, AcAc and BOH levels are determined using an enzymatic assay based on 3-hydroxybutyrate dehydrogenase catalyzed reactions, in which the change in NADH absorbance measured at 340 nm is directly related to either the AcAc and BOH concentrations. This kit can be used various sample types, including serum, plasma, urine, and other biological samples.



Components

This kit is sufficient for up to 100 assays of 3-hydroxybutyric acid (BOH) and up to 100 assays of acetoacetic acid (AcAc) in 96 well plates.

| | |
|------------------------|-------|
| AcAc Buffer | 20 mL |
| Catalog Number MAK134A | |
| BOH Buffer | 20 mL |
| Catalog Number MAK134B | |
| AcAc Reagent | 1 vL |
| Catalog Number MAK134C | |
| BOH Reagent | 1 mL |
| Catalog Number MAK134D | |

AcAc Standard, 80 mM 200 μL
Catalog Number MAK134E

BOH Standard, 80 mM 200 μL
Catalog Number MAK134F

HBDH Enzyme 120 μL
Catalog Number MAK134G

Reagents and Equipment Required but Not Provided.

- Clear 96 well flat-bottom plate suitable for use in UV absorbance assays (Catalog No. CLS3635 or equivalent).
- Spectrophotometric multiwell plate reader

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Briefly centrifuge vials before opening. Use ultrapure water for the preparation of reagents and standards.

AcAc Reagent – Reconstitute with 1,000 μL of water to create a 10 mM solution. Mix well by pipetting, then aliquot and store at -20°C . Use within three weeks of reconstitution.

Storage/Stability

The kit is shipped on dry ice and storage at -20°C , protected from light, is recommended.

Procedures

Serum and plasma samples should be non-hemolyzed and, for best results, assayed immediately. If not assayed, samples can be stored at -80°C for up to 30 days.

Bring all reagents except for the HBDH enzyme to room temperature prior to assay. The HBDH should be kept on ice during use.

AcAc Assay

Standards for AcAc Detection

Add 5 μL of the 80 mM AcAc Standard to 45 μL of water to prepare an 8 mM AcAc Standard solution. Transfer 5 μL of the 8 mM standard and 5 μL of water into separate wells of 96 well plate.

Sample Preparation

Samples can be assayed directly. Aliquot 5 μL of sample into two separate wells of a 96 well plate. One well will be used for the sample activity and one for the sample blank.

AcAc Assay Reaction

1. Set up the AcAc Reaction Mixes according to the scheme in Table 1. 195 μL of the appropriate Reaction Mix is required for each reaction (well).

Table 1.
AcAc Reaction Mixes

| Reagent | Sample and Standards | Sample blank |
|-------------------|----------------------|---------------------|
| AcAc Assay Buffer | 195 μL | 195.5 μL |
| AcAc Reagent | 8 μL | 8 μL |
| HBDH Enzyme | 0.5 μL | – |

2. Add 195 μL of the appropriate Reaction Mix to each of the wells. Mix well using a horizontal shaker or by pipetting, and incubate the reaction at room temperature for 5 minutes. Protect the plate from light during the incubation.
3. Measure the absorbance of the samples and standards at 340 nm (A_{340}).

Results

Calculate the acetoacetic acid (AcAc) concentration from the A_{340} values of the water blank, 8 mM Standard, Sample, and Sample Blank. The concentration of AcAc (mM) in a sample is calculated as:

$$\text{AcAc (mM)} = \frac{A_{\text{sample blank}} - A_{\text{sample}}}{A_{\text{water}} - A_{\text{standard}}} \times 8$$

Note: If the calculated AcAc for the sample is higher than 8 mM, dilute sample in water and repeat the assay. Multiply the results by the dilution factor.

BOH Assay

Standards for BOH Detection

Add 5 μL of the 80 mM BOH Standard to 45 μL of water to prepare an 8 mM BOH Standard solution. Transfer 5 μL of the 8 mM BOH standard and 5 μL of water into separate wells of 96 well plate.

Sample Preparation

Samples can be assayed directly. Aliquot 5 μL of sample into two separate wells of a 96 well plate. One well will be used for the sample activity and one for the sample blank.

BOH Assay Reaction

1. Set up the Reaction Mix according to the scheme in Table 2. 195 μL of the appropriate Reaction Mix is required for each reaction (well).

Table 2.
BOH Reaction Mixes

| Reagent | Sample and Standards | Sample blank |
|------------------|----------------------|---------------------|
| BOH Assay Buffer | 195 μL | 195.5 μL |
| BOH Reagent | 8 μL | 8 μL |
| HBDH Enzyme | 0.5 μL | – |

2. Add 195 μL of the appropriate Reaction Mix to each of the wells. Mix well using a horizontal shaker or by pipetting, and incubate the reaction at room temperature for 15 minutes. Protect the plate from light during the incubation.
3. Measure the absorbance of the samples and standards at 340 nm (A_{340}).

Results

Calculate the 3-hydroxybutyric acid (BOH) concentration from the A_{340} values of the water blank, 8 mM Standard, Sample, and Sample Blank. The concentration of BOH in a sample is calculated as:

$$\text{BOH (mM)} = \frac{A_{\text{sample}} - A_{\text{sample blank}}}{A_{\text{standard}} - A_{\text{water}}} \times 8$$

Note: If the calculated BOH for the sample is higher than 8 mM, dilute sample in water and repeat the assay. Multiply the results by the dilution factor.

Total ketone body (TKB) concentration is calculated as:

$$\text{TKB} = [\text{AcAc}] + [\text{BOH}]$$

Troubleshooting Guide

| Problem | Possible Cause | Suggested Solution |
|--|--|--|
| Assay not working | Omission of step in procedure | Refer and follow Technical Bulletin precisely |
| | Plate reader at incorrect wavelength | Check filter settings of instrument |
| | Type of 96 well plate used | For colorimetric assays, use clear plates |
| Samples with erratic readings | Samples prepared in different buffer | Use the Assay Buffer provided or refer to Technical Bulletin for instructions |
| | Samples used after multiple freeze-thaw cycles | Aliquot and freeze samples if needed to use multiple times |
| | Presence of interfering substance in the sample | If possible, dilute sample further |
| | Use of old or inappropriately stored samples | Use fresh samples and store correctly until use |
| Lower/higher readings in samples and standards | Improperly thawed components | Thaw all components completely and mix gently before use |
| | Use of expired kit or improperly stored reagents | Check the expiration date and store the components appropriately |
| | Allowing the reagents to sit for extended times on ice | Prepare fresh Master Reaction Mix before each use |
| | Incorrect incubation times or temperatures | Refer to Technical Bulletin and verify correct incubation times and temperatures |
| | Incorrect volumes used | Use calibrated pipettes and aliquot correctly |
| Unanticipated results | Samples measured at incorrect wavelength | Check the equipment and filter settings |
| | Samples contain interfering substances | If possible, dilute sample further |
| | Sample readings above the standard | Dilute samples so readings are in range |

KVG,LS,MF,MAM 03/16-1