

Technical Bulletin

Alpha Galactosidase (α -Gal) Activity Assay Kit (Fluorometric)

Catalog Number MAK390

Product Description

Alpha-Galactosidase (α -Gal) hydrolyzes α -galactosyl moieties found in glycolipids and glycoproteins. In mammals, α -Gal hydrolyzes poly- and oligosaccharides commonly found in dietary sources that are difficult to digest. Therefore, α -Gal is used in dietary supplements that help to reduce the production of intestinal gases due to consumption of certain foods. It is known total α -Gal activity is due to two major isozymes with unique, yet different thermostability profiles.

Alpha-Galactosidase A is thermolabile and represents approximately 90% of total α -Gal activity. Fabry Disease, a lysosomal disease disorder, is characterized by mutations in alpha-Galactosidase A. These mutations cause abnormal accumulation of glycosphingolipids in lysosomes.

The Alpha Galactosidase (α -Gal) Activity Assay Kit provides a simple, rapid way to monitor total α -Gal activity in a wide variety of biological samples. In this kit, α -Gal cleaves a synthetic specific substrate, releasing a fluorophore which can be easily quantified ($\lambda_{\text{Ex}} = 360 \text{ nm}/\lambda_{\text{Em}} = 445 \text{ nm}$). The assay is specific (β -galactosidase activity is not detected), sensitive, and capable of detecting as little as $0.1 \mu\text{U}$ of α -Galactosidase activity.

The kit is suitable for the detection of α -Galactosidase activity in tissue homogenates (kidney, etc.), cell lysates (U937, etc.) and biological fluids (saliva, etc.).

Components

The kit is sufficient for 100 fluorometric assays in 96-well plates.

- | | |
|--|-------------------|
| • α -Gal Assay Buffer
Catalog Number MAK390A | 25 mL |
| • α -Gal Stop Buffer
Catalog Number MAK390B | 25 mL |
| • α -Gal Substrate
Catalog Number MAK390C | 220 μL |
| • 4-Methylumbelliferone Standard
Catalog Number MAK390D | 35 μL |
| • α -Gal Positive Control
Catalog Number MAK390E | 1 vial |

Reagents and Equipment Required but Not Provided

- Pipetting devices and accessories (e.g., multichannel pipettor)
- Fluorescence multiwell plate reader
- Flat bottom, white or clear (white is preferred for this assay) 96-well plate. Cell culture or tissue culture treated plates are **not** recommended.
- Refrigerated microcentrifuge capable of $\text{RCF} \geq 12,000 \times g$
- Dounce tissue grinder set (Catalog Number D9063 or equivalent)

MERCK

Precautions and Disclaimer

For Research Use Only. Not for use in diagnostic procedures. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

The kit is shipped on wet ice. Store components at -20 °C, protected from light. Upon opening, use kit within 2 months.

Preparation Instructions

Briefly centrifuge small vials prior to opening.

α -Gal Assay Buffer and Stop Buffer: Store at 2-8 °C or -20 °C. Warm to 37 °C prior to use. Chill an appropriate amount of α -Gal Assay Buffer for use in Sample Preparation

α -Gal Substrate: Light sensitive. Thaw at room temperature. Store at -20 °C.

4-Methylumbelliferone Standard (5 mM): Light sensitive. Thaw at room temperature. Store at -20 °C.

α -Gal Positive Control: Reconstitute with 20 μ L of α -Gal Assay Buffer and mix thoroughly. Store at -20 °C. Keep on ice while in use. Use within two months.

Procedure

Sample Preparation

Tissue and cells

1. Homogenize tissue (10 mg) or pelleted cells ($\sim 5 \times 10^5$ cells) with 100 μ L of ice-cold α -Gal Assay Buffer using a Dounce homogenizer and keep on ice for 10 minutes.
2. Centrifuge samples at $12,000 \times g$ at 4 °C for 10 minutes and collect the supernatant.
3. Dilute the supernatant 10-20-fold in α -Gal Assay Buffer.

4. Add 2-10 μ L of diluted samples into a 96-well plate that will be designated as Sample(s). It is recommended to use 3-5 different volumes of each sample in separate wells to ensure the readings are within the standard curve range and the progress curve rates are within the linear range.
5. Adjust the total volume in each well to 40 μ L with α -Gal Assay Buffer.

Biological Fluids

1. Undiluted fluids can be added directly to the well. Add 2-10 μ L of samples into well(s) in a 96-well plate that will be designated as Samples. It is recommended to use 3-5 different volumes of each sample in separate wells to ensure the readings are within the standard curve range and the progress curve rates are within the linear range.
2. Adjust the total volume in each well to 40 μ L with α -Gal Assay Buffer.

Reagent Background Control

In parallel well(s), prepare Reagent Background Control by adding 40 μ L/well of α -Gal Assay Buffer.

Positive Control

1. Dilute reconstituted α -Gal Positive Control 1:10 with α -Gal Assay Buffer prior to the assay and add 2-6 μ L of diluted α -Gal Positive Control into desired wells(s).
2. Adjust the total volume to 40 μ L/well with α -Gal Assay Buffer.

Note: Do not store unused diluted α -Gal Positive Control.

Standard Curve Preparation

Note: Standards can be prepared at the end of the incubation time (Step 2 of Assay Procedure) and measured in end-point mode.

1. Prepare a 100 μ M 4-Methylumbelliferone (4-MU) Standard by adding 10 μ L of 5 mM 4-MU to 490 μ L of α -Gal Assay Buffer in an amber tube.



2. Prepare a 20 μM 4-MU Standard by diluting the 100 μM Standard solution from Step 1 1:5 by adding 20 μL of 100 μM 4-MU to 80 μL of α -Gal Assay Buffer.
3. Prepare 4-MU Standards according to Table 1. Mix well.

Table 1.
Preparation of 4-MU Standards

Well	20 μM 4-MU Standard	α -Gal Assay Buffer	4-MU (pmol/well)
1	0 μL	50 μL	0
2	2 μL	58 μL	40
3	4 μL	56 μL	80
4	6 μL	54 μL	120
5	8 μL	52 μL	160
6	10 μL	50 μL	200

α -Gal Substrate Working Solution

For each test Sample(s), Positive Control and Reagent Background Control well, prepare 20 μL of α -Gal Substrate Working Solution according to Table 2. Mix well.

Table 2.
Preparation of α -Gal Substrate Working Solution

Reagent	Volume
α -Gal Substrate	2 μL
α -Gal Assay Buffer (pre-warmed to 37 °C)	18 μL

Assay Procedure

1. Add 20 μL of α -Gal Substrate Working Solution to each well containing the test Sample(s), Positive Control, and Reagent Background Control. The total volume in each well (i.e. Samples, Positive Control and Reagent Background Control) should be 60 μL . **Do not add α -Gal Substrate Working Solution to Standard wells.**
2. Mix well and incubate at 37 °C for 2 hours, protected from light.

3. After incubation, add 200 μL of α -Gal Stop Buffer to each well containing Sample(s), Positive Control, Reagent Background Control, and Standards. Mix well.

Measurement

Measure fluorescence intensity (RFU) ($\lambda_{\text{Ex}} = 360 \text{ nm}/\lambda_{\text{Em}} = 445 \text{ nm}$) at 37 °C using end-point mode.

Results

1. Subtract 0 Standard reading (RFU) from all Standard readings.
2. Plot the 4-MU Standard Curve.
3. Subtract the Reagent Background Control reading (RFU) from all Sample readings (RFU).
4. Apply the corrected Sample RFU to the 4-MU Standard Curve to obtain the corresponding pmol of product formed (B). Calculate the activity of α -Galactosidase activity in the sample using the following equation:

$$\alpha\text{-Galactosidase Activity (pmol/h/mg)} \\ \equiv 0.0167 \mu\text{U/mg} =$$

$$\frac{B}{2 \times V \times P} \times D$$

where:

B = 4-MU amount in sample well from Standard Curve (pmol)

2 = Reaction time (hours)

V = Volume of the sample used in the reaction (in mL)

P = Initial Sample Concentration (in mg protein/mL)

D = Sample Dilution Factor (if applicable)

Note: 1 pmol/h = 0.0167 pmol/min
= 0.0167 μU



Unit Definition: One unit of α -Galactosidase activity is the amount of enzyme that generates 1.0 μmol of 4-Methylumbelliferone per minute at pH 4.5 at 37 °C.

Figure 1.
Typical 4-Methylumbelliferone
Standard Curve

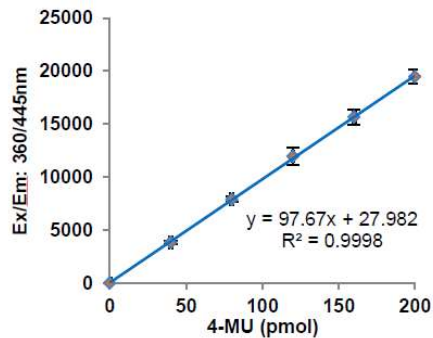


Figure 3.
 α -Galactosidase Activity in undiluted Human
Pooled Saliva (5 μL) measured following the
kit protocol.

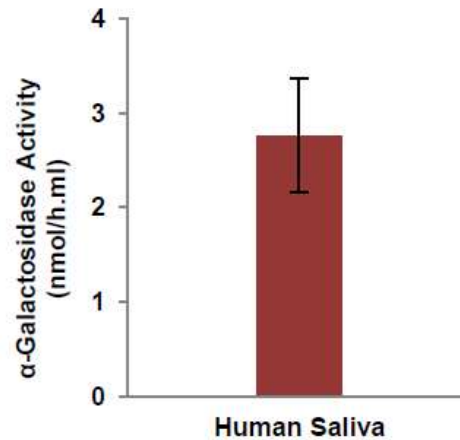
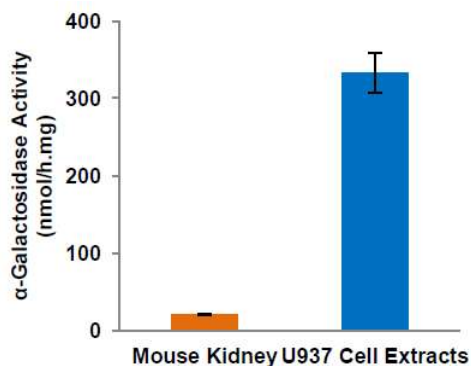


Figure 2.
 α -Galactosidase Activity in Mouse Kidney
Tissue Extracts (1 μg protein) and U937 Cell
Lysates (0.2 μg protein) measured following
the kit protocol.



Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/offices.

Technical Service

Visit the tech service page on our web site at SigmaAldrich.com/techservice.

Standard Warranty

The applicable warranty for the products listed in this publication may be found at SigmaAldrich.com/terms.