

QuantiFluo™ Diamine Oxidase Assay Kit (QFDO-100)

Quantitative Fluorimetric Diamine Oxidase Activity Determination

DESCRIPTION

DIAMINE OXIDASE (DAO) also known as histaminase or amine oxidase (copper containing), is an enzyme involved in the metabolism, oxidation, and inactivation of histamine in animals. Highest levels are observed in the digestive tract and placenta. An imbalance between histamine intake and the capacity for histamine degradation can lead to histamine intolerance (HIT). Measuring DAO activity in serum can be useful in diagnosing HIT.

BioAssay Systems' non-radioactive, fluorimetric DAO assay is based on the oxidation of putrescine to pyrroline plus NH_3 and H_2O_2 . The generated H_2O_2 is then used by HRP to oxidize a dye making it fluorescent. The increase in fluorescence at $\lambda_{\text{ex/em}} = 530/585 \text{ nm}$ is directly proportional to the enzyme activity.

KEY FEATURES

Fast and sensitive. Use of 10 μL sample. Linear detection range 0.5 to 6 U/L for 30 min reaction at 25°C.

Convenient. The procedure involves adding a single working reagent, and reading the fluorescence at 0 and 30 minutes. Room temperature assay. No 37°C heater is needed.

High-throughput. Can be readily automated as a high-throughput 96-well plate assay for thousands of samples per day.

APPLICATIONS

Direct Assays: DAO activity in serum or plasma samples.

KIT CONTENTS (100 TESTS IN 96-WELL PLATES)

Assay Buffer: 10 mL **HRP Enzyme:** 120 μL

Substrate: 120 μL **H_2O_2 Standard:** 100 μL

Dye Reagent: 120 μL

Storage conditions. The kit is shipped at ambient temperature. Store all components at -20°C upon receiving. Shelf life: 6 months after receipt.

Precautions: reagents are for research use only. Briefly centrifuge tubes before opening. Equilibrate all components to room temperature prior to running the assay. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

PROCEDURES

Procedure using 96-well plate

- Internal Standard.** First prepare 500 μL of 8.82 mM H_2O_2 by mixing 5 μL of the H_2O_2 Standard (882 mM) and 495 μL dH_2O . Next mix 20 μL of the 8.82 mM H_2O_2 with 960 μL dH_2O to make a 180 μM internal standard. Use diluted H_2O_2 within 1 hour.
- Prepare sufficient Working Reagent (WR) for all Sample wells by mixing, for each well: 85 μL Assay Buffer, 1 μL HRP Enzyme, 1 μL Substrate and 1 μL Dye Reagent. Prepare sufficient Blank Working Reagent (BWR) for all Sample Blank and Internal Standard wells by mixing for each well, 85 μL Assay Buffer, 1 μL HRP Enzyme, and 1 μL Dye Reagent (i.e. no Substrate).
- Transfer 10 μL of each sample into three separate wells of a black, flat-bottom 96-well plate: one well for Sample measurement (F_S), one for Sample Blank (F_{SB}) and one for the Internal Standard (F_{IS}).
Transfer 10 μL dH_2O to the Sample and Sample Blank wells. Transfer 10 μL of the 180 μM H_2O_2 to the Internal Standard wells.
- Transfer 80 μL WR to each Sample well. Transfer 80 μL BWR to each Sample Blank and Internal Standard well.
- Read fluorescence at $\lambda_{\text{ex/em}} = 530/585 \text{ nm}$ at time 0 and again at time 30 min.

CALCULATION

Subtract the time 0 fluorescence from the time 30 fluorescence for the Sample and Sample Blank wells to compute ΔF_S and ΔF_{SB} respectively. The DAO activity can then be computed as follows:

$$\text{DAO Activity} = \frac{\Delta F_S - \Delta F_{SB}}{F_{IS30} - F_{SB30}} \times \frac{180 \mu\text{M}}{t \text{ (min)}} \times n \text{ (U/L)}$$

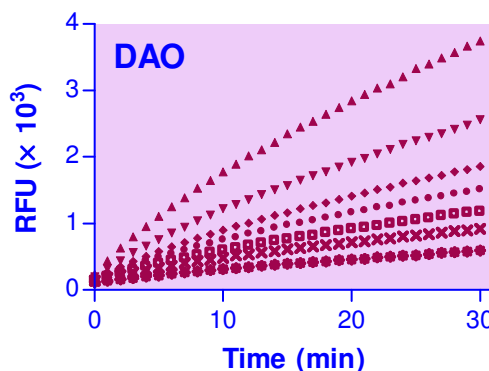
where F_{IS30} and F_{SB30} are the fluorescence readings taken at 30 min for the Internal Standard and Sample Blank respectively and t is the reaction time (30 minutes). n is the sample dilution factor.

Note: if the sample activity is higher than 6 U/L, dilute sample in water and repeat the assay. Multiply the results by the dilution factor. Alternatively, the reaction can be run for a shorter length of time.

Unit definition: 1 Unit (U) of DAO will catalyze the conversion of 1 μmole of putrescine to pyrroline plus NH_3 and H_2O_2 per min at pH 7.5.

MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices, centrifuge tubes, black flat-bottom 96-well plates (e.g. Greiner cat# 655209), and plate reader.



DAO Titration in Human Serum

LITERATURE

- Music, E., et al (2013). Serum diamine oxidase activity as a diagnostic test for histamine intolerance. *Wien Klin Wochenschr* 125:239-243.
- Mondovi B, et al (2013). Effects of amine oxidases in allergic and histamine-mediated conditions. *Recent Pat Inflamm Allergy Drug Discov.* 7(1):20-34.
- Maintz L, Novak N. (2007). Histamine and histamine intolerance. *Am J Clin Nutr.* 85(5):1185-96.

