

Amplite® Colorimetric Tyrosine Assay Kit

Catalog number: 11331
Unit size: 100 Tests

Component	Storage	Amount (Cat No. 11331)
Component A: Tyrosine Enzyme	Freeze (< -15 °C), Minimize light exposure	2 Vials (Lyophilized powder)
Component B: Tyrosine Assay Buffer	Freeze (< -15 °C), Minimize light exposure	1 Bottle (20 mL)
Component C: Tyrosine Standard	Freeze (< -15 °C), Minimize light exposure	1 Vial (Lyophilized powder)

OVERVIEW

Amplite® Colorimetric Tyrosine Assay Kit provides a sensitive and efficient method for quantifying tyrosine in biological and biochemical samples. The assay relies on a colorimetric reaction that generates a measurable chromogenic product proportional to tyrosine concentration, enabling reproducible analysis in research areas such as amino acid metabolism, enzymology, and nutritional science. This assay follows a mix and read format and delivers results in under 60 minutes making it highly convenient for researchers. It supports high-throughput screening and is validated for sample types including plasma, serum, and tissue lysates. The kit includes a tyrosine calibration standard for precise quantification and minimizes interference from common biological substances, ensuring consistent performance across experimental conditions.

AT A GLANCE

Protocol Summary

1. Prepare and add standards and samples (50 µL)
2. Prepare and add Tyrosine working solution to the standards and samples wells (50 µL)
3. Incubate the plate at room temperature for 30 to 60 minutes
4. Monitor the absorbance at 480 nm

Important: Bring all kit components to room temperature before starting the experiment.

KEY PARAMETERS

Absorbance microplate reader

Absorbance 480 nm
Recommended plate Clear plate or White plate/Clear bottom

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

Tyrosine standard solution (100 mM)

1. Add 100 µL Assay Buffer (Component B) to Tyrosine Standard (Component C) and mix well to prepare a 100 mM tyrosine standard solution.

Note: Store any unused Tyrosine stock solution at -20 °C.

Tyrosine Enzyme Stock Solution (50X)

1. Add 100 µL ddH₂O to Tyrosine Enzyme (Component A) and mix

well to prepare a 50X Tyrosine Enzyme stock solutions

Note: Store any unused Tyrosine Enzyme stock solution at -20 °C.

PREPARATION OF STANDARD SOLUTIONS

For convenience, use the Serial Dilution Planner:
<https://www.aatbio.com/tools/serial-dilution/11331>

Preparation of Standard Curve

Add 10 µL of 100 mM Tyrosine Standard solution to 990 µL Tyrosine Assay Buffer (Component B) to prepare a 1 mM Tyrosine solution (Tyr7). Take 300 µL of the Tyr7 solution and perform 1:2 serial dilutions in the Tyrosine Assay Buffer (Component B) to prepare the remaining Tyrosine Standards (Tyr6 to Tyr1), resulting in a concentration range from 500 µM to 15.6 µM.

PREPARATION OF WORKING SOLUTION

Tyrosine working solution

1. Make a 1:50 dilution by adding 20 µL Tyrosine Enzyme Stock Solution (50X) to 1 mL of Tyrosine Assay Buffer (Component B) and mix well.

Note: Tyrosine working solution should be prepared immediately before use. We recommend using the working solution within 2 hours.

SAMPLE EXPERIMENTAL PROTOCOL

Table 1. Layout of Tyrosine standards and test samples in a clear bottom 96-wells microplate. Tyrosine standards (Tyr7-Tyr1 = 1000 µM to 15.6 µM), TS = Test Samples, BL = 0 µM Tyr.

Tyr7	Tyr7	TS	TS				
Tyr6	Tyr6						
Tyr5	Tyr5						
Tyr4	Tyr4						
Tyr3	Tyr3						
Tyr2	Tyr2						
Tyr1	Tyr1						
BL	BL						

The following protocol can be used as a guideline and should be optimized accordingly.

1. Prepare the standards and test samples according to the recommended protocol. Add 50 µL of each into the wells of a microplate.

2. Add 50 μ L Tyrosine working solution to each well containing standards and samples.
3. Incubate the reaction at room temperature for 30 to 60 minutes.
4. Monitor the absorbance with an absorbance plate reader at 480 nm.

EXAMPLE DATA ANALYSIS AND FIGURES

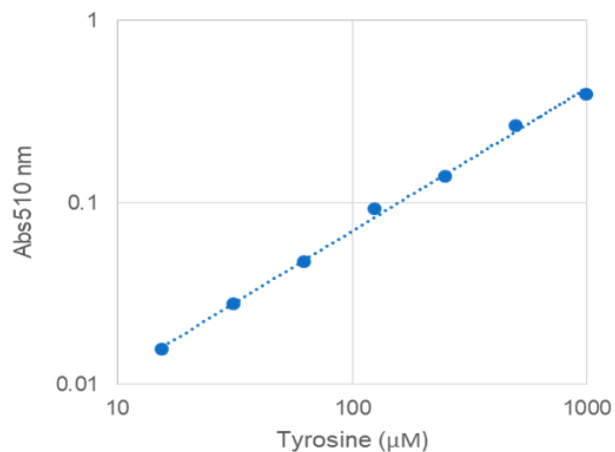


Figure 1. Tyrosine dose response was measured with Amplite® Colorimetric Tyrosine Assay Kit in a 96-well white plate using a SpectraMax microplate reader (Molecular Devices). The signal was acquired at 480 nm.

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 if you have any questions.