

Amplite™ Fluorimetric Formaldehyde Quantitation Kit

Green Fluorescence

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 10057 (200 assays)	Keep at -20 °C, avoid light	Fluorescence microplate readers

Introduction

Formaldehyde is one of the well-identified volatile chemical contaminants responsible for indoor pollution and “building sick” syndrome disease and was recently classified as carcinogenic. The main sources for the HCHO pollution in air include painting, coating material and cigarette smoking. Rapid and accurate measurement of formaldehyde is an important task for biological research, food industry, chemical research and environmental pollution surveillance. There are a few reagents or assay kits available for quantifying formaldehyde. Most of the existing aldehyde test methods are based on separations either by the tedious and expensive HPLC-MS or GC-MS. Our Amplite™ Fluorimetric Formaldehyde Quantitation kit uses a proprietary fluorogenic dye that generates a green fluorescent product upon reacting with formaldehyde. The kit provides a sensitive, one-step fluorometric method to detect as little as 1 μ M of formaldehyde in a 100 μ L assay volume (Figure 1). The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step. Its signal can be easily read by a fluorescence microplate reader at Ex/Em = 410/525 nm.

Kit Key Features

Broad Application:	Quantify formaldehyde in a variety of applications such as enzyme reactions.
Sensitive:	Detect as low as 1 μ M of formaldehyde in a 100 μ L assay volume.
Continuous:	Easily adapted to automation without a separation step.
Convenient:	Formulated to have minimal hands-on time.
Non-Radioactive:	No special requirements for waste treatment.

Kit Components

Components	Amount
Component A: AldeLight™ Green	1 vial
Component B: Assay Buffer	1 bottle (30 mL)
Component C: 37.2% Formaldehyde Standard (12.3 M)	1 vial (100 μ L)
Component D: DMSO	1 vial (100 μ L)

Assay Protocol for One 96-Well Plate

Brief Summary

**Prepare enzyme reaction solution (50 μ L) → Add AldeLight™ Green reaction mixture (50 μ L)
→ Incubate at RT for 20 to 60 minutes → Read Fluorescence at Ex/Em = 410/525 nm**

Note: Thaw all the kit components to room temperature before starting your experiment.

1. Prepare 500X AldeLight™ Green stock solution:

Add 20 μ L of DMSO (Component D) into the AldeLight™ Green vial (Component A) to make 500X stock solution.

Note: The unused AldeLight™ Green solution should be aliquoted, and stored at -20°C (avoid light).

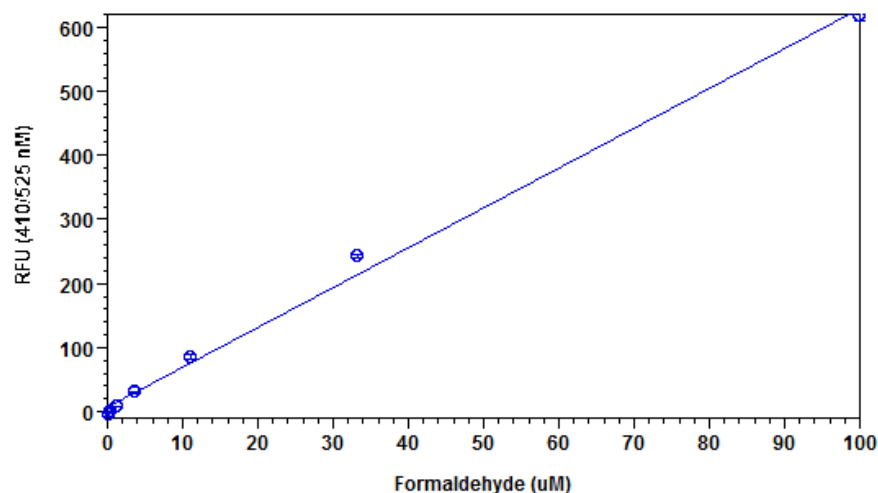


Figure 1. Formaldehyde dose response was measured in a 96-well black plate with Amplitude™ Fluorimetric Formaldehyde Quantitation Kit using a Gemini fluorescence microplate reader (Molecular Devices). As low as 1 μ M of formaldehyde can be detected with 30 minutes incubation time (n=3). *Note: The fluorescence background increases with time, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.*

References

1. Trevor M. Kitson. (1985) High concentrations of aldehydes slow the reaction of cytoplasmic aldehyde dehydrogenase with thiol-group modifiers *Biochem. J.* 228, 765.
2. Crabb DW, Matsumoto M, Chang D, You M (2004). Overview of the role of alcohol dehydrogenase and aldehyde dehydrogenase and their variants in the genesis of alcohol-related pathology. *The Proceedings of the Nutrition Society* 63 (1): 49.
3. Steinmetz CG, Xie P, Weiner H, Hurley TD (1997). Structure of mitochondrial aldehyde dehydrogenase: the genetic component of ethanol aversion. *Structure* 5 (5): 701.
4. O'Donnell JM, Kudej RK, LaNoue KF, Vatner SF, Lewandowski ED. (2004) Limited transfer of cytosolic NADH into mitochondria at high cardiac workload. *Am J Physiol Heart Circ Physiol*, 286, H2237.
5. Zurek G, and karst U (2000). 2,4-Dinitro-3,5,6-trideuterophenylhydrazones for the quantitation of aldehydes and ketones in air samples by liquid chromatography-mass spectrometry. *J of chromatography A*, 869, 251.
6. Ou Z, Ogamo A, Guo L, Konda Y, Harigaya Y, and Nakagawa Y. (1995). Identification and quantitation of choline glycerophospholipids that contain aldehyde residues by fluometric high-performance liquid chromatography. *Analytical biochemistry* 227, 289.

Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.