

Amplite™ Choline Quantitation Kit

Red Fluorescence

Ordering Information:	Storage Conditions:	Instrument Platform:
Product Number: 40007 (200 assays)	Keep in freezer Avoid light	Fluorescence microplate readers

Introduction

Choline is an essential nutrient. Choline and its metabolites play an important role in the structural integrity and signaling of cell membranes and cholinergic neurotransmission (choline synthesis). It is a major source of methyl group via its metabolite, trimethylglycine that participates in the S-adenosylmethionine synthesis pathways. Choline deficiency may cause liver disease, atherosclerosis and possibly neurological disorders. Despite its importance in the central nervous system as a precursor for acetylcholine and membrane phosphatidylcholine, the role of choline in mental illness has been little studied.

Our Amplite™ Choline Quantitation Kit provides one of the most sensitive methods for quantifying choline. The kit uses Amplite™ Red to quantify the concentration of choline, which is related to the production of hydrogen peroxide in the choline oxidase-mediated enzyme coupling reactions. The amount of choline is proportional to the concentration of hydrogen peroxide formed in the enzyme coupling reaction cycle. The kit is an optimized “mix and read” assay that is compatible with HTS liquid handling instruments. It detects as little as 10 picomole choline in 100 µL assay volume (100 nM) as shown in 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 with a fluorescence microplate reader at Ex/Em = ~540/590 nm. Alternatively the assay can also be read at ~576±5 nm with an absorbance microplate reader.

Kit Key Features

Broad Application:	Can be used for quantifying choline in solutions and in cell extracts.
Sensitive:	Detect as low as 10 picomole of choline in solution.
Continuous:	Easily adapted to automation without a separation step.
Convenient:	Formulated to have minimal hands-on time.

Kit Components

Components	Amount
Component A: Amplite™ Red	1 vial
Component B: Choline Probe	2 bottles (lyophilized powder)
Component C: Choline Standard	1 vial (2.8 mg)
Component D: Assay Buffer	1 bottle (25 mL)
Component E: DMSO	1 vial (100µL)

Assay Protocol for One 96-well Plate

Brief Summary

**Prepare choline assay mixture (50 µL) → Add choline standards or choline test samples (50 µL) →
Incubate at RT for 15-60 min → Read fluorescence intensity at Ex/Em = 540/590 nm**

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

Data Analysis

The fluorescence in blank wells (with the assay buffer only) is used as a control, and is subtracted from the values for those wells with the choline reactions. A choline standard curve is shown in Figure 1. The fluorescence background increases with time, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.

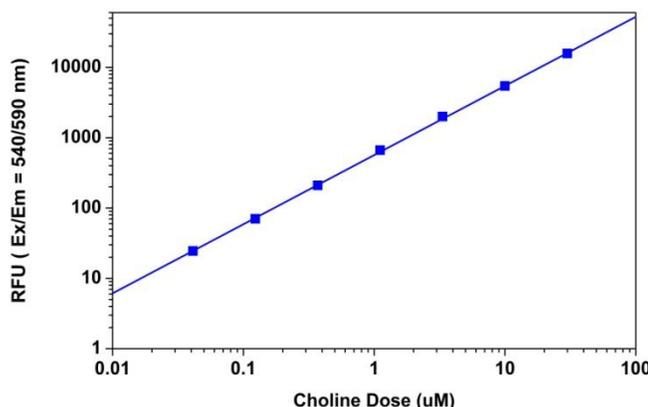


Figure 1. Choline dose response was obtained with Amplite™ Choline Quantitation Kit in a 96-well solid black plate using a Gemini fluorescence microplate reader (Molecular Devices). As low as 100 nM (10 picomole/well) of choline can be detected with 30 minutes incubation time (n=3).

References

1. Al-Saeedi F, Smith T, Welch A. (2007) [Methyl-3H]-choline incorporation into MCF-7 cells: correlation with proliferation, choline kinase and phospholipase D assay. *Anticancer Res*, 27, 901.
2. Park HD, Park KU, Kim KW, Song J, Chang HE, Heo SR, Lee HJ, Kim JQ. (2007) Real-time multiplex PCR assay for genotyping of three apolipoprotein E alleles and two choline acetyltransferase alleles with three hybridization probes. *Clin Chem Lab Med*, 45, 346.
3. Adamczyk M, Brashear RJ, Mattingly PG. (2006) Rapid high-throughput detection of peroxide with an acridinium-9-carboxamide: a homogeneous chemiluminescent assay for plasma choline. *Bioorg Med Chem Lett*, 16, 2407.
4. Shiba K, Ogawa K, Kinuya S, Yajima K, Mori H. (2006) A simple and rapid radiochemical choline acetyltransferase (ChAT) assay screening test. *J Neurosci Methods*, 157, 98.
5. Panfili G, Manzi P, Compagnone D, Scarciglia L, Palleschi G. (2000) Rapid assay of choline in foods using microwave hydrolysis and a choline biosensor. *J Agric Food Chem*, 48, 3403.

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.