

Product datasheet

Flavin Adenine Dinucleotide (FAD) Assay Kit ab204710

1 References 2 Images

Overview

Product name	Flavin Adenine Dinucleotide (FAD) Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Serum, Plasma, Cell Lysate, Cell culture media, Tissue Lysate
Assay type	Enzyme activity
Sensitivity	< 1 nmol/well
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Flavin Adenine Dinucleotide (FAD) Assay Kit (ab204710) is an assay where FAD functions as the cofactor of an oxidase which catalyze the formation of a product that reacts with OxiRed probe generating color and fluorescence. FAD can be detected by either colorimetric (OD=570 nm) or fluorometric (Ex/Em=535/587 nm) methods. The kit provides a rapid, simple, ultra-sensitive, and reliable test suitable for high throughput assay of FAD. The lower limit of detection is less than 1 nM FAD.</p> <p>FAD assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reaction mix - analyze with microplate reader until absorbance reaches 1.8 at 570 nm (or stop before then)
Notes	<p>Flavin Adenine Dinucleotide (FAD) is a redox cofactor which plays an important role in metabolism. FAD exists in different redox states and cycles between FAD, FADH and FADH₂. The primary sources of reduced FAD in eukaryotic metabolism are the citric acid cycle and the beta oxidation reaction pathways.</p>
Platform	Microplate reader

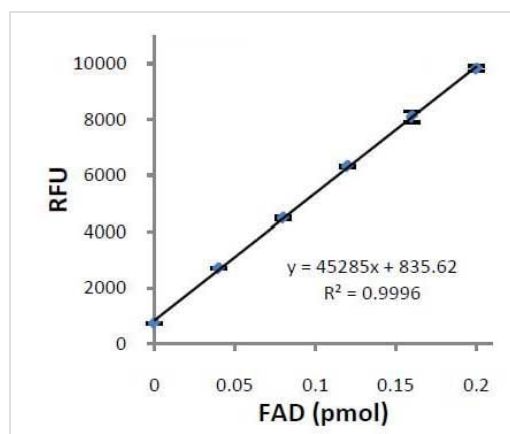
Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
FAD Assay Buffer	1 x 25ml
FAD Enzyme Mix (lyophilized)	1 vial

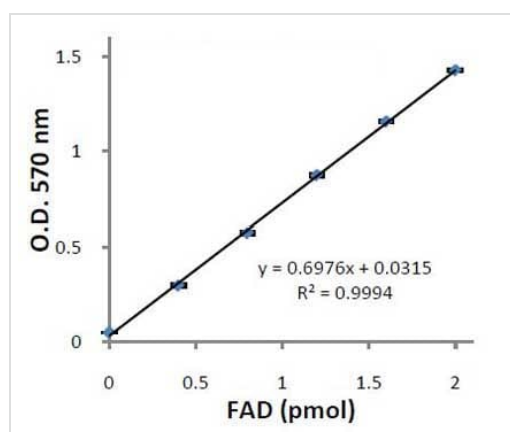
Components	100 tests
FAD Standard (1 nmol, lyophilized)	1 vial
OxiRed Probe (in DMSO)	1 x 200µl
Stop Solution	1 x 1.2ml

Images



Typical FAD Standard Curve (Fluorometric Assay).

Flavin Adenine Dinucleotide (FAD) Assay Kit
(ab204710)



Typical FAD Standard Curve (Colorimetric Assay).

Flavin Adenine Dinucleotide (FAD) Assay Kit
(ab204710)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet

- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors