

Indole Assay Kit ab272536

2 Images

Overview

Product name	Indole Assay Kit
Detection method	Colorimetric
Sample type	Other biological fluids
Assay type	Quantitative
Range	3 μ M - 100 μ M
Product overview	Indole assay kit (272536) is based on a modified version of Ehrlich's and Kovac's reagents, which reacts with indole to produce a colored compound at 565 nm. The intensity of this colored compound is directly proportional to the indole in the sample. Fast and sensitive: Use of 100 μ L sample. Linear detection range from 3 to 100 μ M indole in 96-well plate assay. Convenient: The procedure involves adding a single working reagent, and reading the absorbance immediately.
Tested applications	Suitable for: Functional Studies
Platform	Microplate

Properties

Storage instructions Store at +4°C. Please refer to protocols.

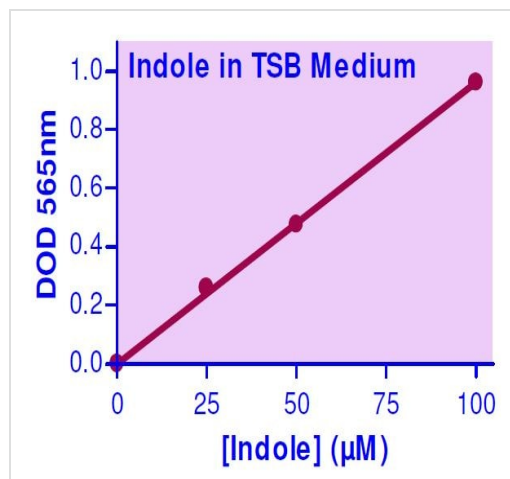
Components	100 tests
Reagent	1 x 12ml
Standard (10 mM Indole)	1 x 100 μ l

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab272536 in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

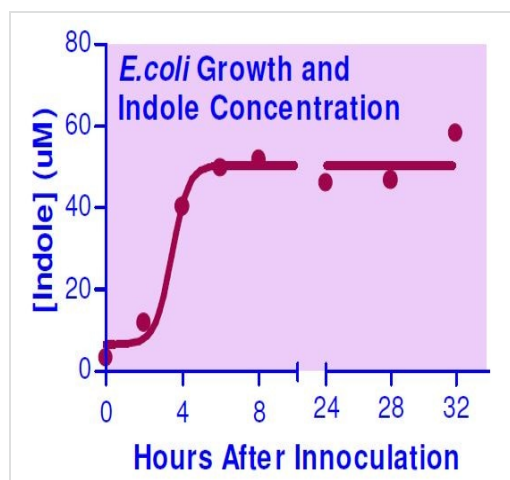
Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration.

Images



Indole Assay Kit Standard Curve

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



E.coli Growth and Indole Concentration

E.coli cells inoculated into 5 mM Tryptophan medium. Medium samples taken every two hours.

Data provided for demonstration purposes only.

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