

AMP Assay Kit (Colorimetric) ab273275

2 Images

Overview

| | |
|------------------|---|
| Product name | AMP Assay Kit (Colorimetric) |
| Detection method | Colorimetric |
| Sample type | Tissue, Adherent cells, Suspension cells |
| Assay type | Quantitative |
| Assay duration | Multiple steps standard assay |
| Product overview | <p>AMP Assay Kit (Colorimetric) (ab273275) provides a convenient method to detect AMP in biological samples. In this assay, AMP is converted to pyruvate in the presence of Pyrophosphate and Phosphoenolpyruvate. This is followed by a set of enzymatic reactions to generate a colored product with a strong absorbance at 570 nm. The absorbance is proportional to the amount of AMP present in samples.</p> <p>The kit is rapid, sensitive, easy to use and high-throughput adaptable. It can measure AMP level lower than 10 µM in various sample types.</p> |
| Notes | <p>This product is manufactured by BioVision, an Abcam company and was previously called K229 AMP Colorimetric Assay Kit. K229-100 is the same size as the 100 test size of ab273275.</p> |
| Platform | Microplate (12 x 8 well strips) |

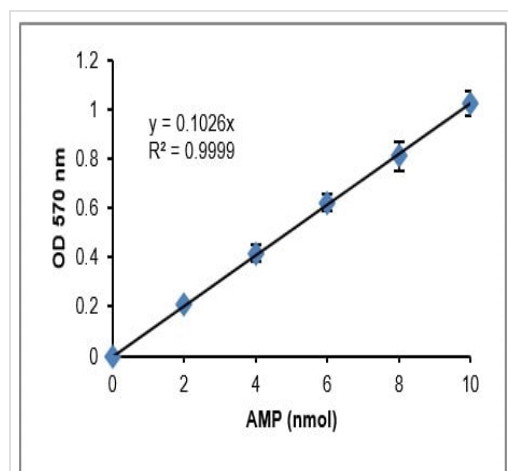
Properties

Storage instructionsStore at -20°C. Please refer to protocols.

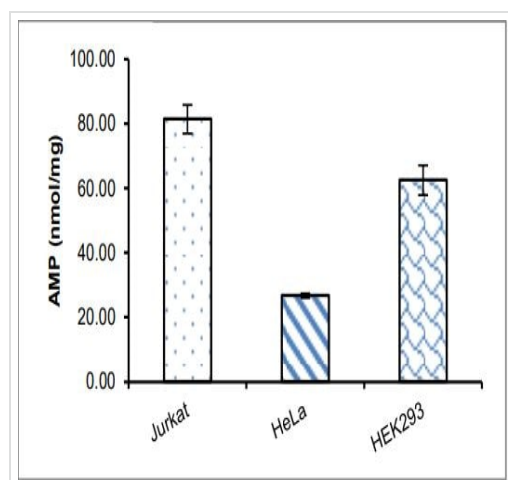
| Components | 100 tests |
|----------------------------|-----------|
| AMP | 1 x 200µl |
| Assay Buffer XV | 1 x 25ml |
| AMP Substrate Mix | 1 vial |
| Developer V | 1 vial |
| Development Enzyme Mix III | 1 x 200µl |
| | |

| Components | 100 tests |
|--------------|-----------|
| OxiRed Probe | 1 x 200µl |

Images



Example data



Example data

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