

Beta Lactamase Activity Assay Kit (Colorimetric) ab197008

[3 Images](#)

Overview

Product name	Beta Lactamase Activity Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Saliva, Urine, Serum, Other biological fluids
Assay type	Enzyme activity
Sensitivity	0.06 mU/well
Assay time	1h 00m
Species reactivity	Reacts with: Mammals, Other species
Product overview	Beta Lactamase Activity Assay Kit (Colorimetric) ab197008 is a simple and sensitive assay that can detect and quantify the enzymatic activity of beta lactamases, a large family of hydrolases.

The beta lactamase assay protocol is based on the hydrolysis of the substrate Nitrocefin, a chromogenic cephalosporin, that results in the generation of a colored product (detectable at OD=490 nm), which is directly proportional to the amount of beta-lactamase activity.

This product can detect as low as 0.06 mU of beta-lactamase activity from bacterial samples or from biological fluids infected with beta-lactamase expressing bacteria.

Beta lactamase assay protocol summary:

- add samples and standards to wells
- add reaction mix
- analyze with a microplate reader in kinetic mode every 2-3 min for 30-60 min

Notes This product is manufactured by BioVision, an Abcam company and was previously called K803 Beta-Lactamase Activity Colorimetric Assay Kit. K803-100 is the same size as the 100 test size of ab197008.

Beta Lactamases (β -Lactamases, β Ls), are a large family of hydrolases comprising more than 850 identified members expressed in Gram-positive and Gram-negative bacteria. β Ls can be classified according to their substrate or inhibitor specificity. These enzymes are capable of hydrolyzing four atom rings known as β -lactams.

Platform

Microplate reader

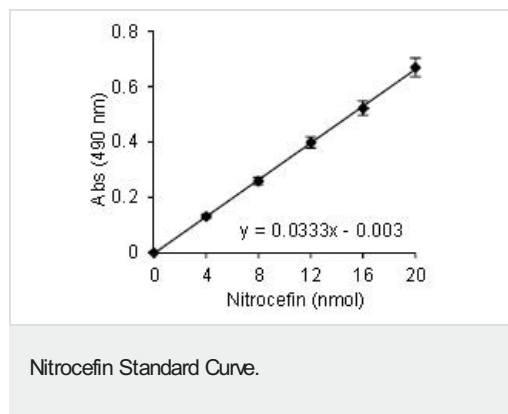
Properties

Storage instructions

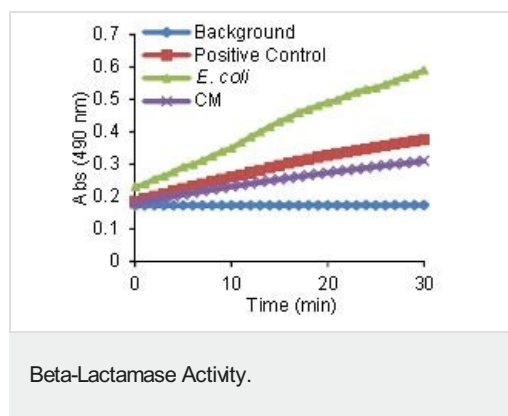
Store at -20°C. Please refer to protocols.

Components	100 tests
Beta Lactamase Assay Buffer	1 x 27ml
Beta Lactamase Hydrolysis Buffer	1 x 100µl
Nitrocefin	1 x 220µl
Positive Control	1 vial

Images

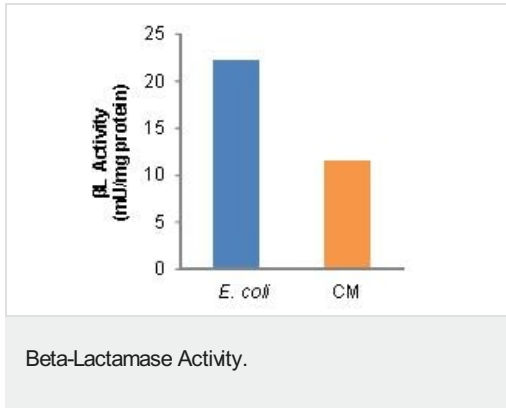


Nitrocefin Standard Curve.



Beta-Lactamase Activity.

Beta-Lactamase activity in *E. coli* culture (5 µl), contaminated media (CM; 30 µl) and Positive Control (4 µl).



Beta-Lactamase Activity of E.coli and contaminated media expressed per milligram of protein.

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