

Product datasheet

Leucine Aminopeptidase (LAP) Activity Assay Kit (Fluorometric) ab234627

[3 Images](#)

Overview

| | |
|-------------------------|---|
| Product name | Leucine Aminopeptidase (LAP) Activity Assay Kit (Fluorometric) |
| Detection method | Fluorescent |
| Sample type | Adherent cells, Suspension cells, Tissue Homogenate |
| Assay type | Enzyme activity |
| Sensitivity | 0.1 mU/well |
| Product overview | Leucine Aminopeptidase (LAP) Activity Assay Kit (Fluorometric) (ab234627) provides a quick, sensitive and easy way for measuring total LAP activity in various samples such as animal tissues or cell cultures. |

In this assay, LAPs hydrolyze leucine from the fluorescent probe and the amount of fluorescent probe detected at Ex/Em 368/460 nm is used to determine the total activity of the LAP enzymes. The assay is simple to perform, high-throughput adaptable and can detect less than 0.1 mU of LAP activity.

Notes This product is manufactured by BioVision, an Abcam company and was previously called K534 Leucine Aminopeptidase (LAP) Activity Assay Kit (Fluorometric). K534-100 is the same size as the 100 test size of ab234627.

Leucine aminopeptidases (EC 3.4.11.1) (LAPs) are a diverse set of exopeptidases that catalyze the hydrolysis of leucine residues from the amino-termini of proteins or peptides. LAPs are ubiquitous enzymes present among animals, plants and prokaryotes. Previously, they were thought to typically play important roles in cell maintenance, growth and development. However, research in the recent years has identified multiple secondary functions for these enzymes in animals and microbes including transcriptional regulation and vesicle transport. Studies have implicated LAP enzymes in tumor cell proliferation, invasion and angiogenesis. Placental LAP is used as a biomarker in ovarian epithelial cancer while adipocyte-derived LAP is used as a marker of endometrial cancer cell proliferation and differentiation.

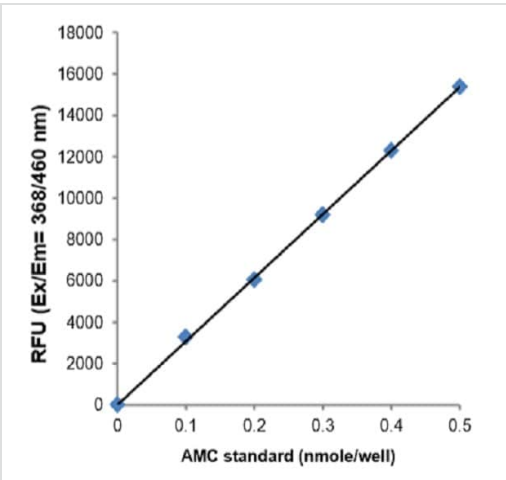
Platform Microplate reader

Properties

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

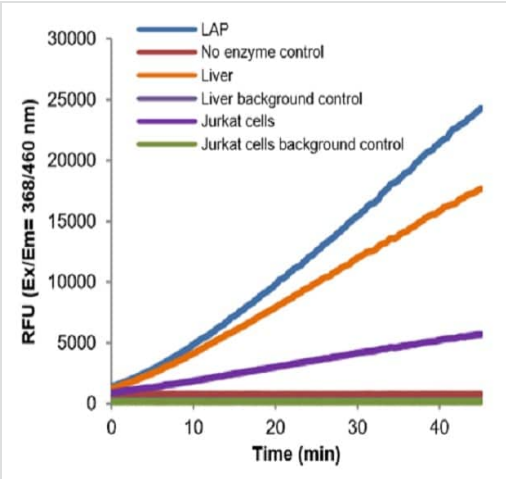
| Components | 100 tests |
|----------------------|-----------|
| AMC Standard | 1 x 100µl |
| LAP Assay Buffer | 1 x 25ml |
| LAP Positive Control | 1 vial |
| LAP Substrate | 1 vial |

Images



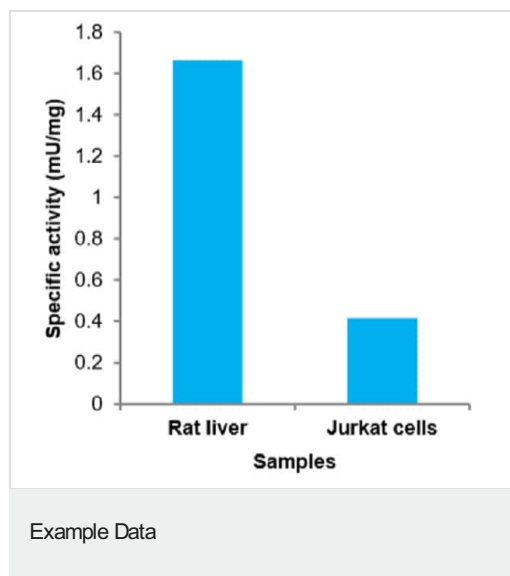
AMC standard curve.

Example data



Reaction kinetics of leucine aminopeptidase activity in rat liver (6.6 µg protein) and Jurkat (human T cell leukemia cell line from peripheral blood) cells (8 µg protein) using appropriate background controls.

Example Data



Leucine aminopeptidase specific activity was calculated in rat liver and Jurkat (Human T cell leukemia cell line from peripheral blood) cell lysates.

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