

Acid Phosphatase Assay Kit (Fluorometric) ab83370

1 图像

概述

产品名称	Acid Phosphatase Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell Lysate, Cell culture media
检测类型	Enzyme activity
灵敏度	0.001 mU/ml
产品概述	Abcam's Acid Phosphatase Assay Kit (Fluorometric) uses non-fluorescent Methylumbelliferyl phosphate disodium (MUP) as the substrate which has Ex/Em=360/440 nm when dephosphorylated by AP. The kit is an ultra-sensitive, simple, direct and HTS-ready fluorometric assay designed to measure AP activity in serum and other bio-samples. Visit our <a href="#">FAQs page</a> for tips and troubleshooting.
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K421 Acid Phosphatase Activity Fluorometric Assay Kit. K421-500 is the same size as the 500 test size of ab83370.</p> <p>The detection sensitivity is ~1 μU, more sensitive than colorimetric assays.</p> <p>The kit is suitable for both research and drug discovery.</p>
平台	Microplate reader

性能

存放说明Store at -20°C. Please refer to protocols.

组件	500 tests
AP Assay buffer	1 x 100ml
AP Enzyme (Lyophilised)	1 vial
MUP Substrate	1 vial
Stop Solution	1 x 25ml

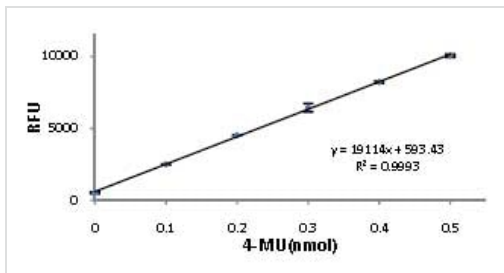
相关性Acid phosphatases (AP) dephosphorylate phosphate groups from phosphate esters under acid conditions. Different acid phosphatase isozymes are found in different organs, and their serum

levels are used as a diagnostic for disease in the corresponding organs. Elevated prostatic acid phosphatase levels may indicate the presence of prostate cancer and elevated tartrate-resistant acid phosphatase levels may indicate bone disease.

## 细胞定位

ACP1: Cytoplasm. ACP2: Lysosome membrane; Single-pass membrane protein. ACP5: Lysosome. ACP6: Isoform 1: Secreted. Isoform 2: Lysosome membrane; Single-pass type I membrane protein.

## 图片



4-Methylumbelliferone (4-MU) Standard Curve.

Functional Studies - Acid Phosphatase Fluorometric Assay Kit (ab83370)

**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.cn/abpromise> or contact our technical team.

## Terms and conditions

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