

ALP Enzyme ab134678

1 References

概述

产品名称	ALP Enzyme
检测类型	Quantitative
产品概述	ALP Enzyme for Alkaline Phosphatase Assay Kit (ab83369) . Reconstitute ALP Enzyme with 1 ml Assay Buffer. <i>The exact concentration/activity of this enzyme is proprietary.</i>

性能

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

组件	标识符	1 units
ALP Enzyme (Lyophilised)	Green	1 vial

相关性 Alkaline phosphatase (ALP) catalyzes the hydrolysis of phosphate esters in alkaline buffer and produces an organic radical and inorganic phosphate. Changes in alkaline phosphatase level and activity are associated with various disease states in the liver and bone.

细胞定位 Cell membrane; Lipid-anchor, GPI-anchor.

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