

# Alkaline Phosphatase Activity Assay Kit II

## Instruction Manual

Catalog Number	PK-CA577-K422															
Description	<p>Alkaline phosphatase (ALP) catalyzes the hydrolysis of phosphate esters in alkaline buffer and produces an organic radical and inorganic phosphate. The change in alkaline phosphatase level and activity is associated with a lot of diseases in the liver and bones. Alkaline phosphatase is also a popular enzyme conjugated to secondary antibody in ELISA.</p> <p>In PromoKine's Alkaline Phosphatase Fluorometric Assay Kit II, ALP cleaves the phosphate group of the non-fluorescent 4-Methylumbelliferyl phosphate disodium salt (MUP) substrate resulting in an intense fluorescent signal (Ex/Em = 360nm/440nm). The kit is an ultra-sensitive, simple, direct and HTS-ready assay designed to measure ALP activity in serum and bio-samples with detection sensitivity ~1 <math>\mu</math>U, more sensitive than colorimetric assays. The kit is suitable for both research and drug discovery.</p>															
Quantity	500 assays															
Applications	<ul style="list-style-type: none"><li>• Measurement of alkaline phosphatase activity</li></ul>															
Sample Type	<ul style="list-style-type: none"><li>• Tissue, cultured cells, media and biological fluids such as serum, plasma and urine</li></ul>															
	<table border="1"><thead><tr><th>Components</th><th>Quantity</th><th>Color Code</th></tr></thead><tbody><tr><td>ALP Assay Buffer</td><td>100 ml</td><td>NM</td></tr><tr><td>MUP Substrate</td><td>1 vial</td><td>Red</td></tr><tr><td>ALP Enzyme</td><td>1 vial</td><td>Green</td></tr><tr><td>Stop Solution</td><td>25 ml</td><td>WM</td></tr></tbody></table>	Components	Quantity	Color Code	ALP Assay Buffer	100 ml	NM	MUP Substrate	1 vial	Red	ALP Enzyme	1 vial	Green	Stop Solution	25 ml	WM
Components	Quantity	Color Code														
ALP Assay Buffer	100 ml	NM														
MUP Substrate	1 vial	Red														
ALP Enzyme	1 vial	Green														
Stop Solution	25 ml	WM														
User Supplied Reagents & Equipment	<ul style="list-style-type: none"><li>• Spectrofluorometer/ Multiwell fluorescence reader</li><li>• 96-well white or black plate with flat bottom</li></ul>															
Storage and Reagents Preparation	<p>Store kit at -20°C, protected from light. Briefly centrifuge all small vials prior to opening. Allow Assay Buffer to warm to room temperature before use. Read entire protocol before performing the assay procedure.</p> <ul style="list-style-type: none"><li>• <b>MUP Solution:</b> Dissolve MUP substrate into 1.2 ml Assay Buffer to generate 5 mM MUP substrate solution. The MUP solution is stable for 2 month at -20°C after dissolved.</li><li>• <b>ALP Enzyme Solution:</b> Reconstitute ALP Enzyme with 1 ml Assay Buffer. The reconstituted enzyme is stable for up to 2 months at 4°C. <b>DO NOT FREEZE!</b></li></ul> <p>Ensure that the Assay Buffer is at room temperature before use. Keep samples and ALP Solution on ice during the assay.</p>															
Assay Protocol	<p><b>1. Sample Preparations:</b> Inhibitors of ALP, like tartrate, fluoride, EDTA, oxalate, and citrate, should be avoided in sample preparation. Serum, plasma, urine, semen, and cell culture media can be assayed directly. Cells (<math>1 \times 10^5</math>) or tissue (~10 mg) can be homogenized in 100 <math>\mu</math>l Assay Buffer, centrifuge to remove insoluble material at 13,000g for 3 minutes. Add test samples directly into 96-well plate, bring total volume to 110 <math>\mu</math>l with Assay Buffer.</p> <p>In order to avoid interference of components in the sample, set a sample background control. Add the same amount of samples into separate wells, bring volume to 110 <math>\mu</math>l. Add 20 <math>\mu</math>l Stop Solution and mix well to terminate ALP activity in the sample.</p> <p><b>2. Reaction Mix:</b> Dilute enough 5 mM MUP substrate solution to 0.5 mM with Assay Buffer (1:10); add 20 <math>\mu</math>l of the 0.5 mM MUP substrate solutions to each well containing the test samples and background controls. Mix well. Incubate the reaction for 30 minutes (or longer if ALP activity in sample is low) at 25°C, protect from light.</p> <p><b>3. Standard Curve:</b> Dilute 10 <math>\mu</math>l of the 5 mM MUP solution with 990 <math>\mu</math>l Assay Buffer to generate 50 <math>\mu</math>M MUP standards. Add 0, 2, 4, 6, 8, 10 <math>\mu</math>l into 96-well plate in duplicate to generate 0, 0.1, 0.2, 0.3, 0.4, 0.5 nmol/well MUP standard. Bring the final volume to 120 <math>\mu</math>l with Assay Buffer.</p> <p>Add 10 <math>\mu</math>l of ALP enzyme solution to each well containing the MUP standard. Mix well.</p> <p>Incubate the reaction for 30 minutes at 25°C, protect from light. The ALP enzyme will convert MUP substrate to equal amount of fluorescent 4-Methylumbelliferone (4-MU).</p> <p><b>4. Measurement:</b> Stop all reactions by adding 20 <math>\mu</math>l Stop Solution into each standard and sample</p>															

reaction except the sample background control reaction (since 20  $\mu$ l Stop Solution has been added into the background control when prepare the sample background control in step 1), gently shake the plate. Measure fluorescence intensity at Ex/Em 360/440 nm using a fluorescence microtiter plate reader.

**5. Calculation:** Correct background by subtracting the value derived from the sample background controls for samples. Plot 4-MU standard Curve. Apply sample readings to the standard curve to get the amount of 4-MU generated by ALP sample. ALP activity of the test samples can be calculated:

**ALP activity= A/V/T (mU/ml)**

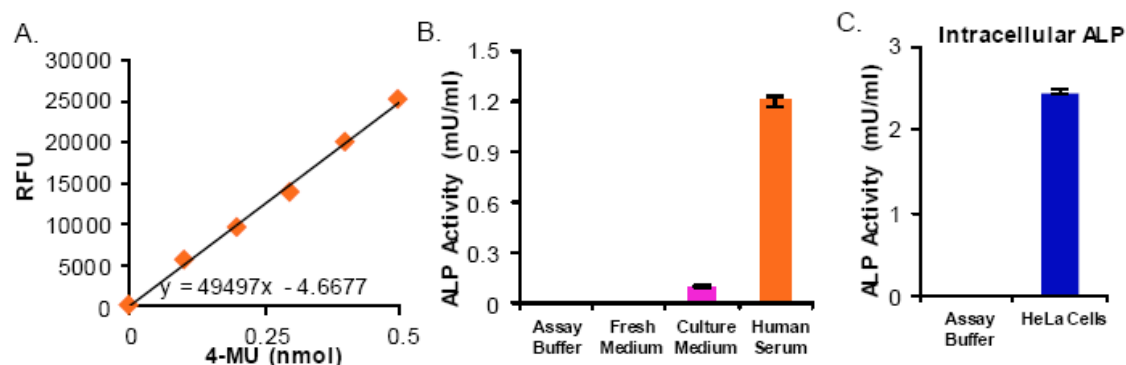
Where: A is amount of 4-MU generated by samples (in nmol).

V is volume of sample added in the assay well (in ml).

T is reaction time (in minutes).

#### Intended Use

For in vitro research use only. Not for diagnostic or therapeutic procedures.



**Figure:** A. 4-MU Standard Curve. B. Measurement of ALP activity in fresh medium (80  $\mu$ l, without culturing), 3-day old HeLa cell cultured medium (80  $\mu$ l) and human serum (80  $\mu$ l, 1:10 diluted). C. Measurement of ALP activity in HeLa cells:  $1 \times 10^4$  HeLa Cells were homogenized, in 200  $\mu$ l of Assay Buffer, diluted 1:10 in Assay Buffer and 80  $\mu$ l was used to measure intracellular ALP activity. Assays were performed according to the kit protocol.

#### PromoCell GmbH

Sickingenstr. 63/65  
69126 Heidelberg  
Germany

Email: [info@promokine.info](mailto:info@promokine.info)  
[www.promokine.info](http://www.promokine.info)

#### USA/Canada

Phone: 1 - 866 - 251 - 2860 (toll free)  
Fax: 1 - 866 - 827 - 9219 (toll free)

#### Deutschland

Telefon: 0800 - 776 66 23 (gebührenfrei)  
Fax: 0800 - 100 83 06 (gebührenfrei)

#### France

Téléphone: 0800 90 93 32 (ligne verte)  
Téléfax: 0800 90 27 36 (ligne verte)

#### United Kingdom

Phone: 0800 - 96 03 33 (toll free)  
Fax: 0800 - 169 85 54 (toll free)

#### Other Countries

Phone: +49 6221 - 649 34 0  
Fax: +49 6221 - 649 34 40