

Product datasheet

# Dihydroxyacetone Phosphate (DHAP) Assay Kit (Fluorometric) ab197003

[2 References](#) [2 Images](#)

Overview

---

<b>Product name</b>	Dihydroxyacetone Phosphate (DHAP) Assay Kit (Fluorometric)
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Serum, Plasma, Other biological fluids, Tissue, Adherent cells, Suspension cells
<b>Assay type</b>	Quantitative
<b>Sensitivity</b>	> 0.5 $\mu$ M
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	Dihydroxyacetone Phosphate (DHAP) Assay Kit (Fluorometric) (ab197003) is suitable for measuring low levels of DHAP typically found in a variety of samples. In this kit, triose phosphate isomerase (TPI) converts dihydroxyacetone phosphate (DHAP) to glyceraldehyde-3-phosphate (GAP), which subsequently undergoes a series of reaction and reduces the probe to generate fluorescence. The fluorescence intensity generated is directly proportional to the amount of dihydroxyacetone Phosphate.

Detection limit: 0.5  $\mu$ M DHAP.

**Notes** This product is manufactured by BioVision, an Abcam company and was previously called K673 PicoProbe™ Dihydroxyacetone Phosphate (DHAP) Fluorometric Assay Kit. K673-100 is the same size as the 100 test size of ab197003.

Dihydroxyacetone Phosphate (DHAP) is an important intermediate in both lipid biosynthesis and glycolysis. In glycolysis, fructose-1,6- diphosphate is converted to dihydroxyacetone phosphate (DHAP) and glyceraldehyde-3-phosphate (GAP) by aldolase. Both DHAP and GAP serve as the intracellular pool for triose phosphate. DHAP can be further converted into GAP by Triose Phosphate Isomerase (TPI).

In humans, TPI deficiency is a rare autosomal disease. It causes hemolytic anemia, neurological diseases, and even death due to blockage of the glycolytic pathway and accumulation of DHAP in erythrocytes.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

**Platform** Microplate reader

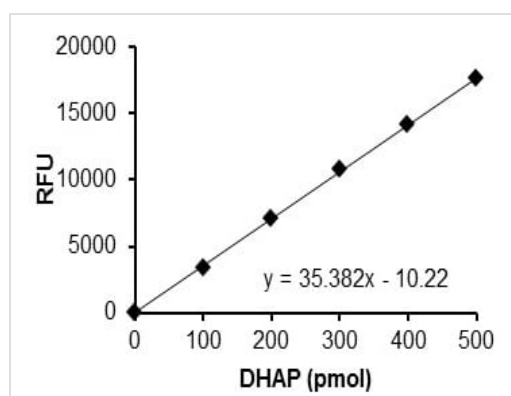
## Properties

### Storage instructions

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

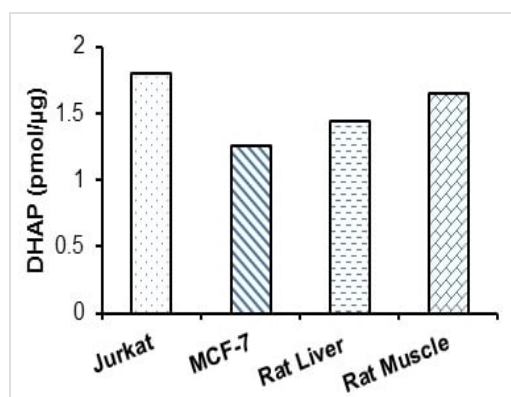
Components	Identifier	100 tests
DHAP Assay Buffer		1 x 25ml
DHAP Developer		1 vial
DHAP Enzyme Mix		1 vial
DHAP Standard		1 vial
PicoProbe	Blue	1 x 400µl

## Images



Typical DHAP Standard Curve obtained following assay protocol.

Functional Studies - Dihydroxyacetone Phosphate  
(DHAP) Assay Kit (Fluorometric) (ab197003)



Measurement of DHAP level in a variety of samples: Jurkat (250 µg), and MCF-7 (150 µg) cell lysate, and in rat liver (50 µg) and rat muscle (150 µg) tissue lysate.

Functional Studies - Dihydroxyacetone Phosphate  
(DHAP) Assay Kit (Fluorometric) (ab197003)

**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