

6-Phosphofructokinase Activity Assay Kit (Colorimetric) ab155898

[22 References](#) [1 Image](#)

Overview

Product name	6-Phosphofructokinase Activity Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Cell Lysate, Tissue Lysate
Assay type	Enzyme activity
Sensitivity	< 1 mU/well
Range	2 nmol/well - 10 nmol/well
Assay time	1h 30m
Species reactivity	Reacts with: Mammals, Other species
Product overview	6-Phosphofructokinase Activity Assay kit ab155898 is simple, robust, and rapid and can detect phosphofructokinase activity less than 1 mU.

In the PFK activity assay protocol, PFK (phosphofructokinase) converts fructose-6-phosphate and ATP to fructose-diphosphate and ADP. The ADP in the presence of substrate and enzyme mix is converted to AMP and NADH, which reduces a colorless probe to a colored product with strong absorbance at 450 nm.

PFK activity assay protocol summary:

- add samples and standards to wells
- add reaction mix
- analyze with a microplate reader over 20-60 min

Notes This product is manufactured by BioVision, an Abcam company and was previously called K776 Phosphofructokinase (PFK) Activity Colorimetric Assay Kit. K776-100 is the same size as the 100 test size of ab155898.

Phosphofructokinase (PFK) is a key glycolytic enzyme and plays a major regulatory role during glycolysis. This enzyme is present in bacteria, plants and animals. There are 2 types of PFKs - PFK1 and PFK2. In the presence of ATP, PFK1 & PFK2 catalyzes the conversion of fructose-6-phosphate to fructose-1,6-diphosphate and fructose-2,6-diphosphate respectively and ADP. PFK has 3 major isoforms in mammals: PFK-M (muscle), PFK-L (liver) and PFK-P (platelet).

Platform Microplate reader

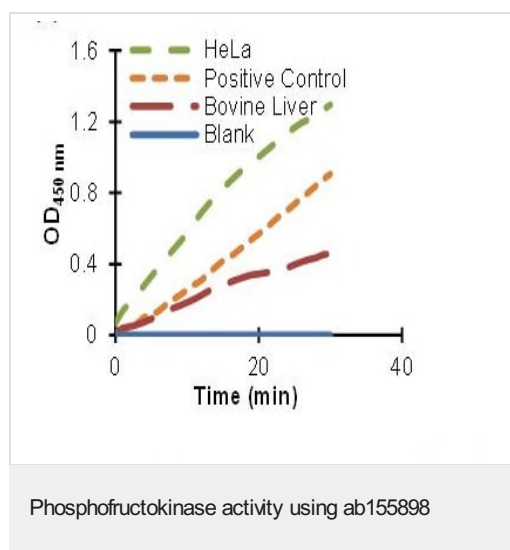
Properties

Storage instructions

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

Components	100 tests
Assay Buffer VII	1 x 27ml
ATP III	1 vial
CK Enzyme Mix	1 vial
Developer Solution III	1 vial
NADH Standard I	1 vial
PFK Positive Control	1 vial
PFK Substrate	1 vial

Images



PFK activity determined in bovine liver and HeLa cell lysate. Assay performed following kit protocol.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Promise 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