

Product datasheet

# 6-Phosphofructokinase Activity Assay Kit (Colorimetric) ab155898

1 Image

Overview

---

<b>Product name</b>	6-Phosphofructokinase Activity Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell Lysate, Tissue Lysate
<b>Assay type</b>	Enzyme activity
<b>Sensitivity</b>	< 1 mU/well
<b>Range</b>	2 nmol/well - 10 nmol/well
<b>Assay time</b>	1h 30m
<b>Species reactivity</b>	<b>Reacts with:</b> Other species, Mammals
<b>Product overview</b>	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** 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

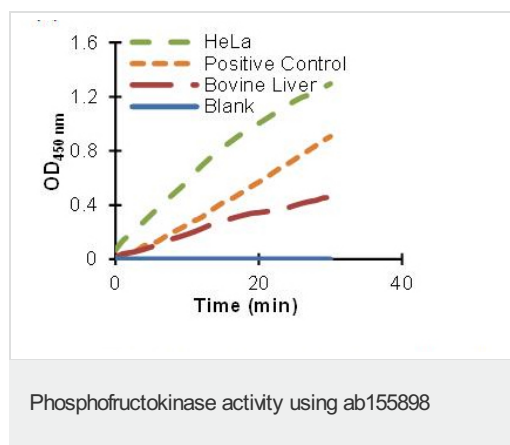
Properties

---

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	Identifier	100 tests
ATP	Orange	1 vial
NADH Standard	Yellow	1 vial
PFK Assay Buffer	WM	1 x 27ml
PFK Developer	Red	1 vial
PFK Enzyme Mix	Green	1 vial
PFK Positive Control	Purple	1 vial
PFK Substrate	Blue	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 AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## 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