

Product datasheet

Gamma Glutamyl Transferase (GGT) Assay Kit (Fluorometric) ab241030

★★★★★ 2 Abreviews 2 Images

Overview

Product name	Gamma Glutamyl Transferase (GGT) Assay Kit (Fluorometric)
Detection method	Fluorescent
Species reactivity	<b>Reacts with:</b> Mammals
Product overview	Gamma Glutamyl Transferase (GGT) Assay Kit (Fluorometric) (ab241030) provides a convenient tool for sensitive detection of the GGT in a variety of samples.

The GGT in sample will recognize L-γ-Glutamyl-AMC as a specific substrate leading to proportional fluorescence development. The activity of GGT can be easily quantified fluorometrically (Ex/Em = 365/460 nm).

This assay detects GGT activity as low as 0.02 mIU in sample.

Notes	This product is manufactured by BioVision, an Abcam company and was previously called K785 Gamma Glutamyl Transferase (GGT) Activity Fluorometric Assay Kit. K785-100 is the same size as the 100 test size of ab241030.
-------	--

Properties

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

Components	100 tests
AMC Standard	1 x 100µl
GGT Assay Buffer	1 x 25ml
GGT Positive Control	1 vial
GGT Substrate	1 unit

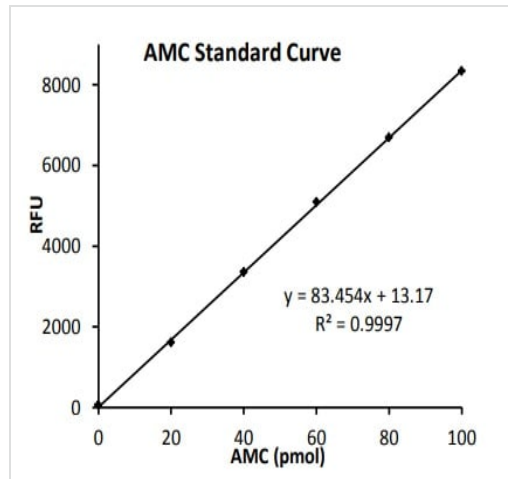
Relevance	Glutamyltransferase (gGT) from human kidney is a membrane bound enzyme that transfers the gamma glutamly moiety of gamma glutamyl compounds such as glutathione to an acceptor, which may be an amino acid or a di or tripeptide or water. The principal functions of gGT may be
-----------	--

hydrolysis and metabolism of glutathione. Various human gGT isoforms have been found. The molecular mass of gGT for normal kidney is 90kDa and gGT can be used as a renal cell marker.

## Cellular localization

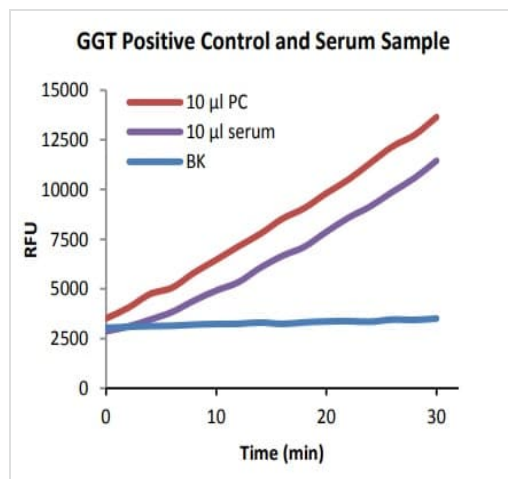
Cell Membrane

## Images



AMC Standard Curve.

AMC Standard Curve.



GGT Positive Control and Human Serum Sample.

GGT Positive Control and Serum Sample.

**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