

# Lipase for Triglyceride Quantification Kit ab89001

## 3 References

### 概述

产品名称	Lipase for Triglyceride Quantification试剂盒
产品概述	ab89001 is the catalytic component of Abcam's <b>Triglyceride Quantification Kit (ab65336)</b> .

### 性能

**存放说明** Store at -20°C. Please refer to protocols.

组件	1 units
Lipase (Lyophilized)	1 vial

**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.cn/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors