

Enzyme Mix for Triglyceride Quantification Kit ab155160

製品の概要

製品名	Enzyme Mix for Triglyceride Quantification Kit
製品の概要	Enzyme mix for Triglyceride Quantification Kit (ab65336)

製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
Enzyme Mix VI	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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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