



Product datasheet

Triglyceride Standard ab103967

概述

产品名称	Triglyceride Standard
产品概述	Kit component from Triglyceride Quantification Kit (ab65336) .
说明	Frozen storage may cause the triglyceride standard to separate from the aqueous phase. To re-dissolve, keep the cap tightly closed, place in a hot water bath (~80-100°C) for 1 minute or until the standard looks cloudy, vortex for 30 seconds, the standard should become clear. Repeat the heat and vortex one more time. The triglyceride standard is now completely in solution, and ready to use.
经测试应用	适用于: Purification

性能

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

组件	300 µl
Triglyceride Standard	1 x 300µl

应用

The Abpromise guarantee [Abpromise™](#)承诺保证使用ab103967于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
Purification		Use at an assay dependent concentration.

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