

## Product datasheet

### Cholesterol Assay Buffer ab136652

#### Overview

<b>Product name</b>	Cholesterol Assay Buffer
<b>Product overview</b>	ab136652 is the Cholesterol Assay Buffer for <b>ab65390</b> HDL and LDL/VLDL Cholesterol Assay Kit

#### Properties

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

Components	25 ml
Cholesterol Assay Buffer	1 x 25ml

**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