

# 2X LDL/VLDL Precipitation Buffer AB105138

## 1 References

### Overview

<b>Product name</b>	2X LDL/VLDL Precipitation Buffer
<b>Assay type</b>	Quantitative
<b>Product overview</b>	2X LDL/VLDL Precipitation Buffer for <a href="#">ab65390 HDL and LDL/VLDL Cholesterol Assay Kit</a> .

### Properties

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

Components	Identifier	10 ml
2X LDL/VLDL Precipitation Buffer	NM	1 x 10ml

**Relevance** HDL (high-density-lipoprotein), LDL (low-density lipoprotein) and VLDL (very low-density lipoprotein) are lipoproteins that act as carrier proteins for cholesterol. The regulation of HDL-cholesterol and LDL-cholesterol plays a central role in various disease development. It is well known that low levels of HDL and high level of LDL are associated with an increased risk of cardiovascular events.

**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