

Product datasheet

10x Red Blood Cell (RBC) Lysis Buffer ab204733

8 References

Overview

Product name	10x Red Blood Cell (RBC) Lysis Buffer
Sample type	Tissue, Whole Blood
Product overview	10x Red Blood Cell (RBC) Lysis Buffer (ab204733) provides a quick and efficient method of lysing red blood cells. Human whole blood is composed of 45% red blood cells. Without the removal of red blood cells, it is difficult to analyze the phenotype and function of leukocytes in whole blood. Haemoglobin and other red blood cell contents can also interfere with several chemical assays.

Notes

This product is manufactured by BioVision, an Abcam company and was previously called 5831 10X Red Blood Cell Lysis Buffer. 5831-100 is the same size as the 100 ml size of ab204733.

10x Red Blood Cell (RBC) Lysis Buffer uses ammonium chloride method to lyse red blood cells without affecting leukocytes, normal tissue, or tumor cells.

Properties

Storage instructions Store at room temperature. Please refer to protocols.

Components	100 ml
10X Red Blood Cell (RBC) Lysis Buffer	1 x 100ml

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