

Product datasheet

Blocking Buffer 10X ab210904

2 References

Overview

Product name Blocking Buffer 10X

Product overview Blocking Buffer 10X is recommended for use with the Matched Antibody Pair kits during ELISA development. The Blocking Buffer 10X is a BSA based buffer (10 %w/v) that is designed to reduce background interference due to non-specific binding.

The 10X Blocking Buffer should be diluted to 1X using 1X PBS before use.

See our [website](#) for a protocol.

The Blocking Buffer 10X is also available as part of our Matched Antibody Pair Kit Accessory Pack ([ab210905](#)).

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	50 ml
Blocking Buffer 10X (ab210904)	1 x 50ml

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