

Product datasheet

ELISA Coating Buffer 1X ab210899

1 References

Overview

Product name ELISA Coating Buffer 1X

Product overview ELISA Coating Buffer 1X is recommended for use with the Matched Antibody Pair kits for plate coating during ELISA development. See our website for more protocol instructions.

ELISA Coating Buffer 1X is supplied ready to use.

The ELISA Coating Buffer 1X 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	120 ml
ELISA Coating Buffer 1X (ab210899)	1 x 120ml

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