

ELISA Coating Buffer 1X ab210899

2 References

概述

产品名称	ELISA Coating缓冲液1X
产品概述	<p>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.</p> <p>ELISA Coating Buffer 1X is supplied ready to use.</p> <p>The ELISA Coating Buffer 1X is also available as part of our Matched Antibody Pair Kit Accessory Pack (ab210905).</p>

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors