

ELISA Accessory Pack ab210905

2 References

概述

产品名称	ELISA Accessory Pack
产品概述	<p>ELISA Accessory Pack is recommended for use with the Matched antibody pair kits for sandwich ELISA.</p> <p>See our website for a protocol.</p> <p>Please store all components at 4°C, apart from the Streptavidin HRP Solution which needs to be stored at -20°C. The plate can be stored at 4°C or at room temperature. You can also refer to the datasheets for the individual components for storage information.</p>

性能

存放说明 Please refer to protocols.

组件	1 packs
10X Wash Buffer PT (ab206977)	1 x 1000ml
96-Well Microplate	10 x 96 tests
Blocking Buffer 10X (ab210904)	1 x 50ml
ELISA Coating Buffer 1X (ab210899)	1 x 120ml
Stop Solution (ab210900)	1 x 120ml
Streptavidin HRP Solution (ab210901) 50 µg/mL	1 x 110µl
TMB Substrate (ab210902)	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