

Antibody Diluent 5BI ab221825

SimpleStep ELISA

概述

产品名称

抗体稀释剂5BI

产品概述

Antibody diluent to be used with some of our SimpleStep® ELISA kits.

Supplied ready to use.

性能

存放说明

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

组件	6 ml
Antibody Diluent 5BI	1 x 6ml

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