

## Product datasheet

### Antibody Diluent 5BI ab221825

SimpleStep ELISA<sup>®</sup>

#### 製品の概要

製品名	Antibody Diluent 5BI
製品の概要	Antibody diluent to be used with some of our SimpleStep <sup>®</sup> ELISA kits.  Supplied ready to use.

#### 製品の特性

保存方法	Store at +4°C. Please refer to protocols.
------	---

内容	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.co.jp/abpromise> or contact our technical team.

#### Terms and conditions

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