

### 10X Wash Buffer PT ab206977

SimpleStep ELISA

#### 製品の概要

---

製品名	10X Wash Buffer PT
製品の概要	Abcam's 10X Wash Buffer PT is a used with our range of SimpleStep ELISAs and Matched Antibody Pair Kits.

#### Instructions for use

10X Wash Buffer PT should be diluted to 1X before use. To prepare 1X Wash Buffer PT dilute 10X Wash Buffer PT with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL 10X Wash Buffer PT with 45 mL deionized water. Mix thoroughly and gently.

#### 製品の特性

---

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

**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