

## Product datasheet

### Sample Diluent NBP ab221826

SimpleStep ELISA

#### Overview

**Product name** Sample Diluent NBP

**Product overview** Sample diluent to be used with some of our SimpleStep<sup>®</sup> ELISA kits.

Supplied ready to use or can be diluted to appropriate concentration.

Note: This product contains bovine plasma (citrate).

#### Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	20 ml
Sample Diluent NBP	1 x 20ml

**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.com/abpromise> or contact our technical team.

## **Terms and conditions**

---

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