

Pre-Coated 96 Well Microplate (12 x 8 well strips) SimpleStep ELISA® ab206978

SimpleStep ELISA

製品の概要

製品名	Pre-Coated 96 Well Microplate (12 x 8 well strips) SimpleStep ELISA®
製品の概要	<p>Abcam's SimpleStep ELISA Pre-Coated 96 Well Microplate (12 x 8 well strips) is a component used with our range of SimpleStep ELISAs.</p> <p>The microplate is pre-coated with the monoclonal antibody of a proprietary tag and is only suitable for use with our SimpleStep ELISA range.</p>

製品の特性

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

内容	1 units
Plate Seals	1 unit
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit

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