

PDH Immunocapture Kit ab109802

2 References 1 图像

概述

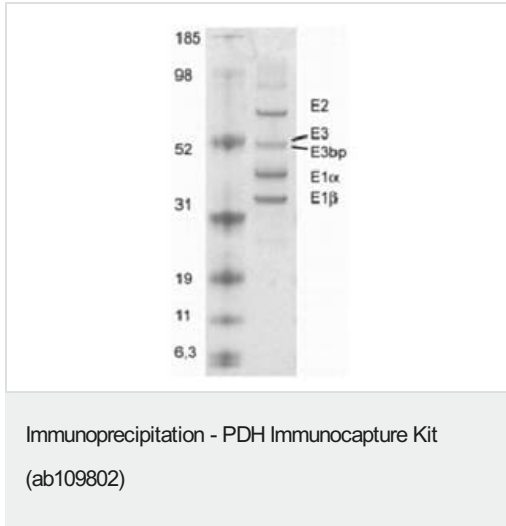
产品名称	PDH免疫捕获试剂盒
检测类型	Quantitative
种属反应性	与反应: Cow, Human
产品概述	250 µg monoclonal antibody irreversibly crosslinked to protein G-agarose beads which can immunocapture PDH complex from heart mitochondria.
Note: The immunocapture protocol for this kit requires Abcam reagent detergent lauryl maltoside ( <a href="#">ab109857</a> /MS910).	

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	250 µg
monoclonal antibody irreversibly crosslinked to protein G-agarose beads	1 x 250µg

图片



PDH immunoprecipitation using antibody **ab109866**/MSP01c crosslinked to protein G-agarose beads as product ab109802

**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