

Complex IV Immunocapture Kit ab109801

11 References **1 图像**

概述

产品名称	Complex IV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	<p>ab109801 Complex IV Immunocapture Kit allows isolation of the cytochrome c oxidase complex from small amounts of tissue.</p> <p>250 µg, 500 µg or 750 µg monoclonal antibodies irreversibly crosslinked to protein G-agarose beads which can immunocapture ~25 µg, ~50 µg or ~75 µg respectively of Complex IV from heart mitochondria.</p> <p>This kit facilitates subsequent analysis of both assembly state and accrued post-translational modifications of all 13 subunits of Complex IV. The immunoprecipitated Complex IV shows cytochrome c oxidase activity while bound to the beads and this activity is fully sensitive to cyanide. Uses for ab109801 include but are not limited to examining alterations of Complex IV subunits in inherited mitochondrial diseases, Alzheimer's disease, schizophrenia and cancer.</p> <p>Note: The immunocapture protocol for this kit requires Abcam detergent lauryl maltoside (ab109857/MS910).</p>

说明

Related products

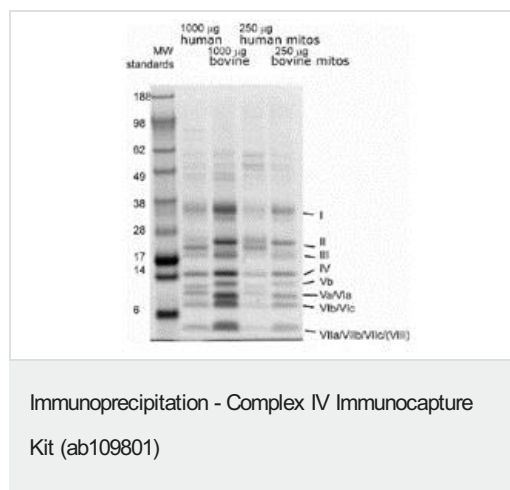
Review the [mitochondrial assay guide](#), or the full [metabolism assay guide](#) to learn about more assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also how to assay metabolic function in live cells using your plate reader.

性能

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

组件	250 µg
Monoclonal Complex IV antibody irreversibly crosslinked to protein G-agarose beads	1 x 250µg

图片



Complex IV immunoprecipitation using antibody

ab109863/MS101c crosslinked to protein G-agarose beads as product ab109801

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