

Complex III Immunocapture Kit ab109800

5 References [1 图像](#)

概述

产品名称	Complex III免疫捕获试剂盒
检测类型	Quantitative
种属反应性	与反应: Mouse, Rat, Cow, Human
产品概述	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 III from heart mitochondria.

ab109800 Complex III Immunocapture Kit allows isolation of the ubiquinol-cytochrome c oxidoreductase complex from small amounts of tissue. This facilitates subsequent analysis of assembly state, activity and the extent of post translational modifications including oxidative damage that occur with aging. Uses for ab109800 include research on genetic mitochondrial disease, where complex III may have an additional role in the stabilization of complex I.

Note: The immunocapture protocol for this kit requires Abcam detergent lauryl maltoside ([ab109857/MS910](#)).

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.

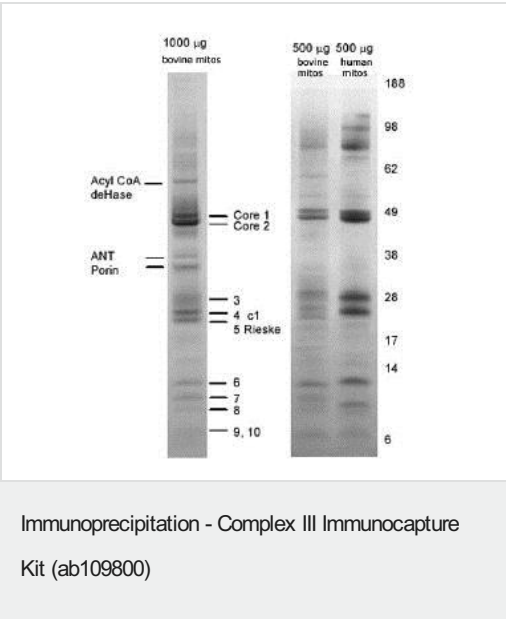
性能

组件	250 µg
monoclonal Complex III antibody irreversibly cross-linked to protein G-agarose beads	1 x 250µg

相关性

Complex II (succinate-ubiquinone oxidoreductase), one of the mitochondrial respiratory chain complexes, transfers electrons from succinate generated during the citric acid cycle to Complex III (ubiquinolcytochrome c oxidoreductase), via a mobile electron shuttle, ubiquinone. Complex III transfers electrons to Complex IV (cytochrome c oxidase) via another mobile electron shuttle, cytochrome c.

图片



Complex III immunoprecipitation using antibody **ab109862**/MS301 crosslinked to protein G-agarose beads as product ab109800.

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