

Complex II Immunocapture Kit ab109799

7 References [1 图像](#)

概述

产品名称	Complex II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 ~13 µg, ~26 µg or ~39 µg respectively of Complex II from heart mitochondria. Isolated mitochondria are available from several species for use as control samples.

The Complex II Immunocapture Kit allows isolation of the succinate dehydrogenase 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 the Complex II immunocapture kit include research on aging genetic mitochondrial disease, and various cancers due to mutations in this enzyme complex such as paragangliomas and phaeo-chromocytomas.

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

存放说明

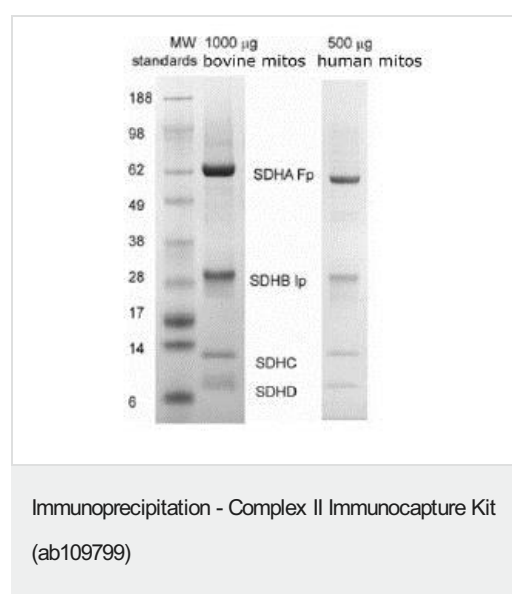
Store at +4°C. Please refer to protocols.

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

相关性

Complex II is also called succinate ubiquinone oxidoreductase or more commonly succinate dehydrogenase complex. This complex is composed of four nuclear encoded subunits and contains a flavin (FAD), non-heme iron centers and a b-type cytochrome as prosthetic groups. It is both a component of the electron transport chain and an enzyme of the Krebs cycle. Complex II deficiencies are seen in OXPHOS genetic disease and found in a type of cancer called paraganglioma.

图片



Complex II immunoprecipitation using antibody **ab109865**

crosslinked to protein G-agarose beads as product ab109799.

Complex II was immunoprecipitated from a lauryl maltoside detergent extract of 1000 ug bovine heart mitochondria or 500 ug human heart mitochondria. The gel was stained with coomassie brilliant blue.

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