

Product datasheet

Complex III Immunocapture Kit ab109800

[3 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	<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 III from heart mitochondria.</p> <p>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.</p> <p>Note: The immunocapture protocol for this kit requires Abcam detergent lauryl maltoside (ab109857/MS910).</p>

经测试应用 **适用于: IP**

性能

组件	250 µg	500 µg	750 µg
monoclonal Complex III antibody irreversibly cross-linked to protein G-agarose beads	1 x 250µg	1 x 500µg	1 x 750µg

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.

应用

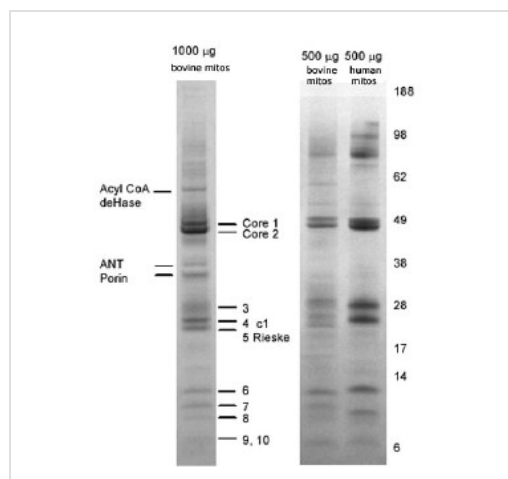
Our [Abpromise guarantee](#) covers the use of **ab109800** in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
----	------	----

IP

Use at an assay dependent dilution.

图片



Complex III immunoprecipitation using antibody [ab109862/MS301](#) crosslinked to protein G-agarose beads as product [ab109800](#).

Immunoprecipitation - Complex III Immunocapture Kit ([ab109800](#))

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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