

ATP Synthase Immunocapture Kit ab109715

[10 References](#) [1 Image](#)

Overview

Product name	ATP Synthase Immunocapture Kit
Assay type	Quantitative
Species reactivity	Reacts with: Rat, Cow, Human
Product overview	250 µg of monoclonal antibodies irreversibly crosslinked to protein G-agarose beads which can immunocapture up to ~125 µg, ~250 µg or ~500 µg respectively of ATP synthase (Complex V) from heart mitochondria.

ab109715 allows isolation of the ATP synthase complex from small amounts of tissue. This facilitates subsequent analysis of assembly state and activity. Thus the enzyme retains oligomycin sensitive ATP hydrolysis activities after isolation. Finally, the extent of post translational modifications including oxidative damage can be readily analyzed by proteomic approaches or antibody detection of these modifications. Potential uses for ab109715 include but are not limited to examining alterations of ATP synthase subunits in inherited mitochondrial diseases, Alzheimer's disease, angiogenesis, hypertension and aging.

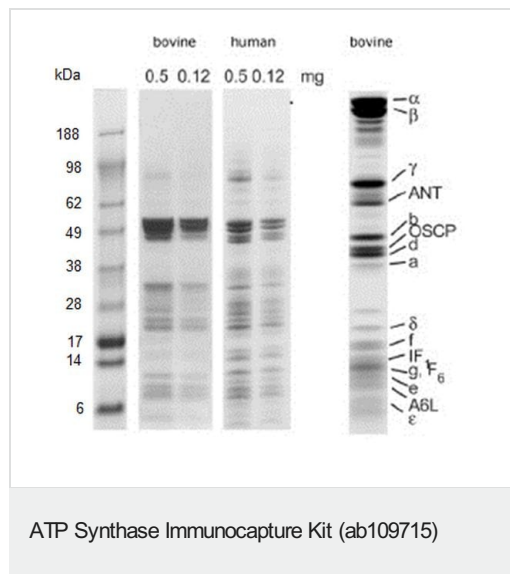
Note: The immunocapture protocol for this kit requires Abcam detergent lauryl maltoside ([ab109857](#)/MS910).

Properties

Storage instructions	Store at +4°C. Please refer to protocols.
-----------------------------	---

Components	250 µg
ATP synthase monoclonal antibody coupled agarose beads	2 x 250µg

Images



Using the immunoprecipitation protocol provided, ATP synthase was isolated from various samples by antibody **ab109867** crosslinked to protein G-agarose beads (ab109715). Protein bands identities were confirmed by mass spectrometry.

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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors