

Adenine Nucleotide Translocase (ANT) Immunocapture Kit ab109713

[5 References](#) [画像数 1](#)

製品の概要

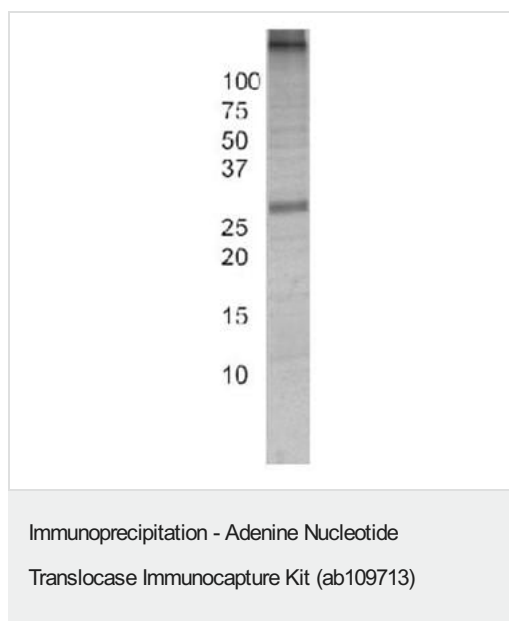
製品名	Adenine Nucleotide Translocase (ANT) Immunocapture Kit
アッセイタイプ	Quantitative
種交差性	交差種: Cow, Human
製品の概要	Capture monoclonal antibody irreversibly crosslinked to protein G-agarose beads which can immunocapture Adenine Nucleotide Translocase from heart mitochondria. Note: The immunocapture protocol for this kit requires Abcam detergent lauryl maltoside (ab109857 /MS910).

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	250 µg
Adenine Nucleotide Translocase monoclonal antibody coupled agarose Beads	1 x 250µg

画像



Adenine Nucleotide Translocase immunoprecipitation using antibody **ab109864**/MS101c crosslinked to protein G-agarose beads as product ab109713

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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