

Product datasheet

Adenine Nucleotide Translocase (ANT)  
Immunocapture Kit ab109713

5 References 1 Image

Overview

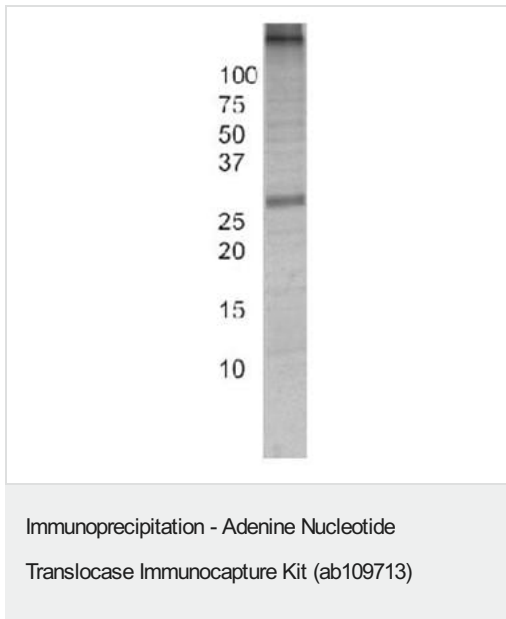
Product name	Adenine Nucleotide Translocase (ANT) Immunocapture Kit
Assay type	Quantitative
Species reactivity	<b>Reacts with:</b> Cow, Human
Product overview	Capture monoclonal antibody irreversibly crosslinked to protein G-agarose beads which can immunocapture Adenine Nucleotide Translocase from heart mitochondria.  Note: The immunocapture protocol for this kit requires Abcam detergent lauryl maltoside ( <a href="#">ab109857</a> /MS910).

Properties

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

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

Images



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.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