

Mitochondrial DNA Isolation Kit ab65321

★★★★★ 2 Abreviews 56 References

Overview

Product name	Mitochondrial DNA Isolation Kit
Sample type	Tissue, Adherent cells, Suspension cells
Assay type	Direct
Assay time	3h 00m
Product overview	<p>Mitochondrial DNA Isolation Kit ab65321 provides convenient tools for isolating mtDNA from a variety of cells and tissues in high yield and purity, without contaminations from genomic DNA. The purified mtDNA can be used for a variety of studies such as enzyme manipulations, Southern blotting, cloning, PCR analysis, and amplifications</p>

Please note: One kit is sufficient for 50 samples if Adherent cell or Suspension cells are used and for 25 samples if tissue is used.

Mitochondrial DNA isolation protocol summary:

- collect cells by centrifugation at 600 x g for 5 min and wash with PBS
- centrifuge for 5 min and resuspend pellet in cytosol extraction buffer
- incubate on ice for 10 min
- homogenize in dounce tissue grinder
- centrifuge for 10 min at 700 x g to pellet nuclei and intact cells
- transfer supernatant and centrifuge at 10,000 g for 30 min
- remove supernatant, resuspend pellet in cytosol extraction buffer and centrifuge at 10,000 g for 30 min
- resuspend pellet (mitochondria) and lyse in mitochondrial lysis buffer for 10 min
- add enzyme mix and incubate for 60 min
- add ethanol and incubate for 10 min
- centrifuge for 5 min and retain pellet (mitochondrial DNA)

Notes This product is manufactured by BioVision, an Abcam company and was previously called K280 Mitochondrial DNA Isolation Kit. K280-50 is the same size as the 1 kit size of ab65321.

Mitochondria are semi-autonomous organelles which functions in aging process, apoptosis, anti-HIV drugs, and cancers. Mitochondrial DNA (mtDNA) has a very high mutation rate and the mutations on mtDNA appear to be related to certain diseases such as diabetes, Alzheimer's disease, and muscle disorders. Isolation and quantification of mtDNA are often required to study the relationships between the diseases and mtDNA.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of

products that contain European Authorisation list (Annex XIV) substances.
It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	1 kit
5X Cytosol Extraction Buffer I	1 x 20ml
Enzyme B Mix	1 vial
Lysis Buffer III	1 x 1.8ml
TE Buffer I	1 x 1.5ml

Relevance Mitochondria are semiautonomous organelles which function in the aging process, apoptosis, anti-HIV drugs, and cancers. Mitochondrial DNA (mtDNA) has a very high mutation rate and the mutations on mtDNA appear to be related to certain diseases such as diabetes, Alzheimer's disease, and muscle disorders.

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