

Product datasheet

Yeast nuclei isolation kit ab206997

[1 Image](#)

Overview

Product name Yeast nuclei isolation kit

Product overview Abcam's Yeast nuclei isolation kit (ab206997) provides an easy and fast method for isolation of nuclei from yeast cells for downstream applications such as DNA-protein interaction, RNA-protein interaction and protein-protein interaction studies, DNase I foot-printing analysis, enzymatic assays, pull-down assays, as well as western blot and ELISA.

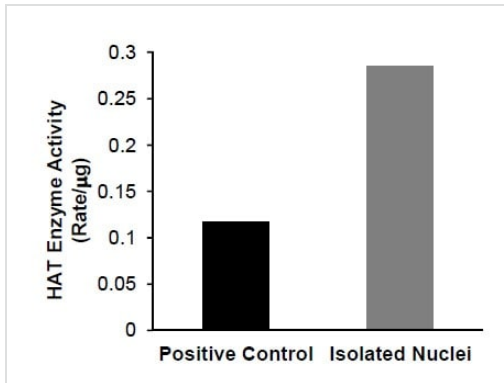
To understand the nature of the replication and expression of the yeast genome, it is essential to have a method for the preparation of nuclei at different stages of growth. Abcam's Yeast nuclei isolation kit enables easy and fast purification of nuclei from yeast cells using yeast cell wall lysis and homogenization.

Properties

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

Components	50 tests
1 M DTT	1 x 1ml
Buffer A	1 x 50ml
Buffer B	1 x 50ml
Buffer C	1 x 90ml
Lysis Enzyme Cocktail	1 x 500µl
Protease Inhibitor Cocktail (Lyophilized) (22mg)	1 vial

Images



Functional Activity of Isolated Yeast Nuclei. Purified yeast nuclei were analyzed for HAT enzyme activity using H3 substrate. In yeast, H3-dependent HAT activity is nuclear specific. Positive Control: Nuclear Extract (NE, 4 mg/mL). Purified yeast nuclei show significant HAT Enzyme Activity. Yeast nuclei were isolated following the kit protocol.

Functional Activity of Isolated Yeast Nuclei.

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