

## Product datasheet

# Preadipocyte Isolation kit ab196988

★★★★★ 1 Abreviews 1 References 2 Images

### Overview

<b>Product name</b>	Preadipocyte Isolation kit
<b>Sample type</b>	Tissue
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse, Rat, Other species, Mammals <b>Predicted to work with:</b> Mammals

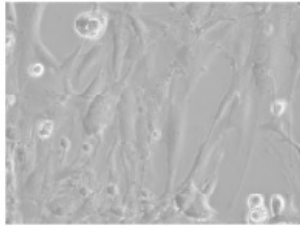
**Product overview** Preadipocyte Isolation kit (ab196988) provides the reagents and tools to isolate the stromal vascular fraction from adipose tissue. The resulting stromal vascular fraction is cultured on a tissue culture plate, to which the preadipocytes adhere. The preadipocytes retain the ability to proliferate and differentiate into adipocytes when treated with differentiation-inducing components. The preadipocytes can be used to study the process of adipogenesis, while the differentiated adipocytes can be used to study lipolysis, endocrine activity, cell-signaling and metabolic dysfunction.

### Properties

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

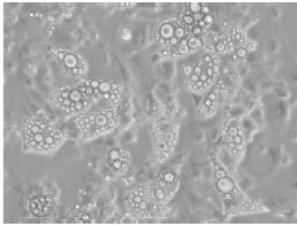
Components	Identifier	5 g
Cell Strainer (100 µm)		10 units
Cell Strainer (70 µm)		10 units
Collagenase	NM, blue	1 x 10ml
Collagenase Stop Buffer	NM	1 x 90ml
PBS		1 x 90ml
Red Blood Cell Lysis Buffer		1 x 10ml

### Images



Example data

Preadipocytes from mouse adipose tissue after 4 days of culturing.



Example data

Adipocytes 7 days after differentiation from preadipocytes using a 3T3-L1 Differentiation Kit. Visible lipid droplets confirm the isolation of preadipocytes.

**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