

# Hematopoietic Stem Cell Marker (CD34, CD59, CD90 / Thy1, CD38, c-Kit) Antibody Panel - Human ab254022

リコンビナント

1 References [画像数 10](#)

## 製品の概要

製品名	Hematopoietic Stem Cell Marker (CD34, CD59, CD90 / Thy1, CD38, c-Kit) Antibody Panel - Human
種交差性	交差種: Human
製品の概要	<p>Hematopoietic Stem Cell Marker (CD34, CD59, CD90 / Thy1, CD38, c-Kit) Antibody Panel - Human ab254022 contains multiple trial-sized versions of anti-human antibody clones against CD34, CD59, CD90 / Thy1, CD38, c-Kit, specifically selected for high performance in various applications. This panel contains 5 recombinant rabbit monoclonal antibodies against human CD34, CD59, CD90 / Thy1, CD38, c-Kit. They are provided as a sampler panel to allow you to easily evaluate each in your required applications.</p> <p>For guidelines on how to use each antibody within the panel, please consult the individual datasheet for each antibody.</p> <p>Panel contains:</p> <ul style="list-style-type: none"> <li>- Rabbit monoclonal [EP373Y] to CD34 (20 µL) <a href="#">ab81289</a></li> <li>- Rabbit monoclonal [EPR6425(2)] to CD59 (20 µL) <a href="#">ab133707</a></li> <li>- Rabbit monoclonal [EPR3133] to CD90 / Thy1 (20 µL) <a href="#">ab133350</a></li> <li>- Rabbit monoclonal [EPR4106] to CD38 (20 µL) <a href="#">ab108403</a></li> <li>- Rabbit monoclonal [YR145] to c-Kit (20 µL) <a href="#">ab32363</a></li> </ul>

**特記事項**

**Explore our range of antibody sample panels** designed to provide you with a variety of trial-size antibodies in a convenient and cost-effective format.

**Carrier-free formulations** of our recombinant antibodies are available and ready to use for

multiplex IHC analysis including Imaging Mass Cytometry™. Please refer to the 'Associated products' section below.

Our RabMAb® technology is a patented hybridoma-based technology for making rabbit monoclonal antibodies. For details on our patents, please refer to [RabMAb® patents](#).

## 製品の特性

### 保存方法

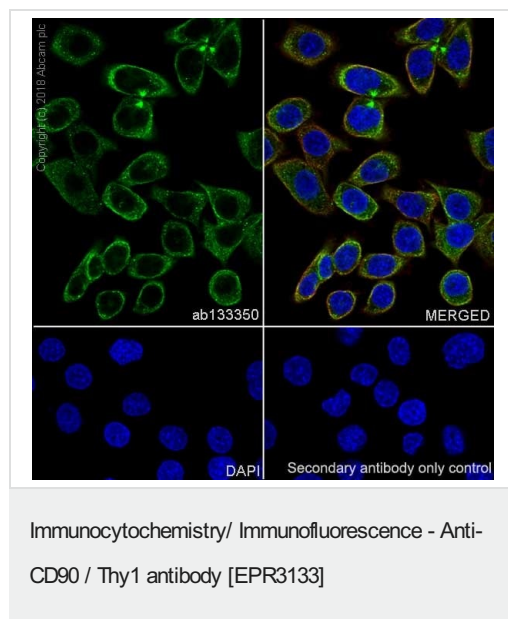
Store at -20°C. Please refer to protocols.

内容	1 kit
<a href="#">ab81289 - Anti-CD34 antibody [EP373Y]</a>	2 x 10µl
<a href="#">ab108403 - Anti-CD38 antibody [EPR4106]</a>	2 x 10µl
<a href="#">ab133707 - Anti-CD59 antibody [EPR6425(2)]</a>	2 x 10µl
<a href="#">ab133350 - Anti-CD90 / Thy1 antibody [EPR3133]</a>	2 x 10µl
<a href="#">ab32363 - Anti-c-Kit antibody [YR145]</a>	2 x 10µl

### 細胞内局在

CD34: Membrane. CD38: Membrane. CD59: Cell membrane. Secreted. Soluble form found in a number of tissues. c-Kit: Cell membrane and Cytoplasm. Detected in the cytoplasm of spermatozoa, especially in the equatorial and subacrosomal region of the sperm head. CD90 / Thy1: Cell membrane.

## 画像

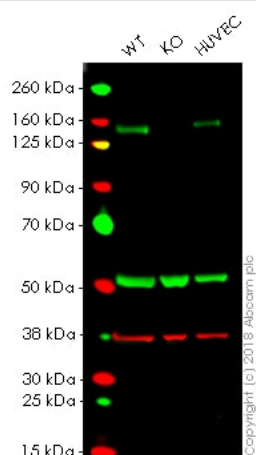


[ab133350](#) staining CD90 / Thy1 in the SH-SY5Y human

neuroblastoma epithelial cell line by ICC/IF

(Immunocytochemistry/immunofluorescence). Cells were fixed with 4% Paraformaldehyde and permeabilized with 0.1% Triton X-100.

Samples were incubated with primary antibody (1/100). [ab150077](#) an Alexa Fluor® 488-conjugated Goat anti-rabbit IgG (1/1000) was used as the secondary antibody. DAPI was used as a nuclear counter stain and Ab195889 Anti-alpha Tubulin antibody [DM1A] - Microtubule Marker (Alexa Fluor® 594) was used to counter stain microtubules.



Western blot - Anti-c-Kit antibody [YR145]

**All lanes :** Anti-c-Kit antibody [YR145] ([ab32363](#)) at 1/1000 dilution

**Lane 1 :** Wild-type HAP1 whole cell lysate

**Lane 2 :** KIT knockout HAP1 whole cell lysate

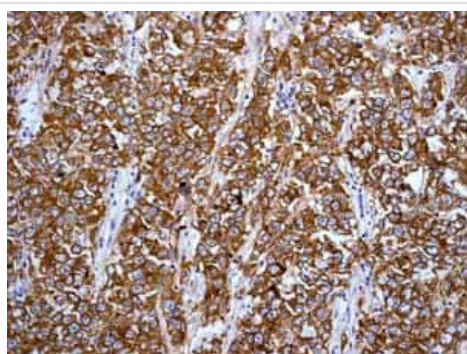
**Lane 3 :** HUVEC whole cell lysate

Lysates/proteins at 40 µg per lane.

**Predicted band size:** 110 kDa  
Lanes 1 - 3: Merged signal (red and green). Green - [ab32363](#) observed at 109 kDa. Red - loading control, [ab9484](#), observed at 37 kDa.

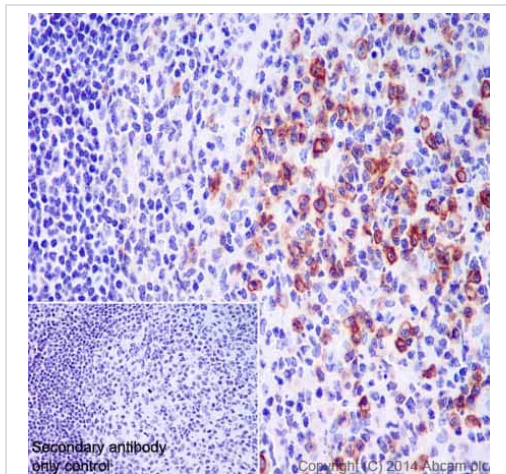
**Lanes 1 - 3:** Merged signal (red and green). Green - [ab32363](#) observed at 109 kDa. Red - loading control, [ab9484](#), observed at 37 kDa.

[ab32363](#) was shown to recognize c-Kit in wild-type HAP1 cells as signal was lost at the expected MW in KIT knockout cells. Additional cross-reactive bands were observed in the wild-type and knockout cells. Wild-type and KIT knockout samples were subjected to SDS-PAGE. Ab32363 and [ab9484](#) (Mouse anti-GAPDH loading control) were incubated overnight at 4°C at 1/1000 dilution and 1/20000 dilution respectively. Blots were developed with Goat anti-Rabbit IgG H&L (IRDye® 800CW) preabsorbed [ab216773](#) and Goat anti-Mouse IgG H&L (IRDye® 680RD) preabsorbed [ab216776](#) secondary antibodies at 1/20000 dilution for 1 hour at room temperature before imaging.



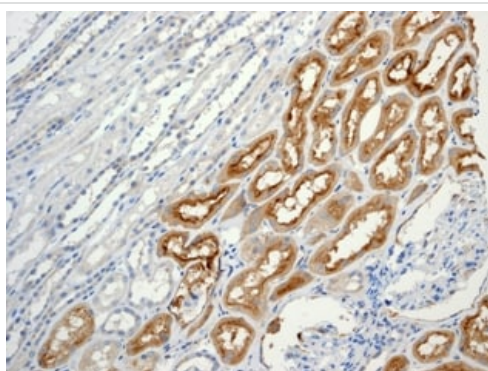
Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti-c-Kit antibody [YR145]

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) analysis of human seminoma tissue labelling c-Kit with unpurified [ab32363](#).



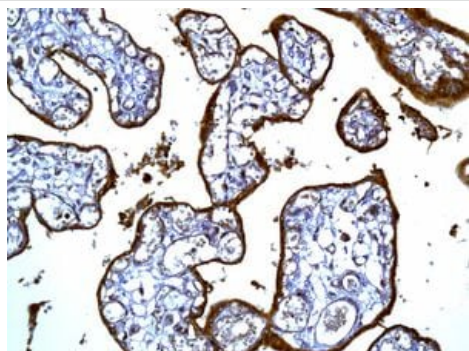
Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti-CD38 antibody  
[EPR4106]

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) analysis of human tonsil tissue labelling CD38 with purified **ab108403** at 1/500. Heat mediated antigen retrieval was performed using Tris/EDTA buffer pH 9. **ab97051**, a HRP-conjugated goat anti-rabbit IgG (H+L) was used as the secondary antibody (1/500). Negative control using PBS instead of primary antibody. Counterstained with hematoxylin.



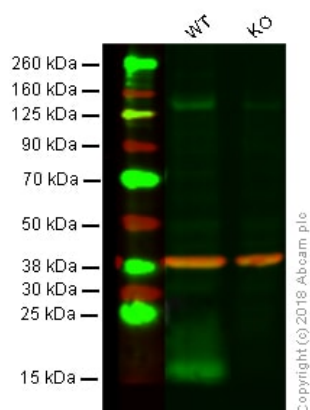
Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti-CD90 / Thy1 antibody  
[EPR3133]

Immunohistochemical analysis of paraffin-embedded Human normal kidney tissue labelling CD90 / Thy1 with **ab133350** at 1/100 dilution.



Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti-CD59 antibody [EPR6425(2)]

Immunohistochemical analysis of CD59 in paraffin embedded Human placenta tissue labelled with **ab133707** at a 1/100 dilution.



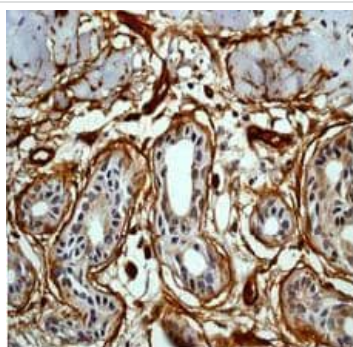
Western blot - Anti-CD59 antibody [EPR6425(2)]

**Lane 1:** Wild-type HAP1 whole cell lysate (40 µg)

**Lane 2:** CD59 knockout HAP1 whole cell lysate (40 µg)

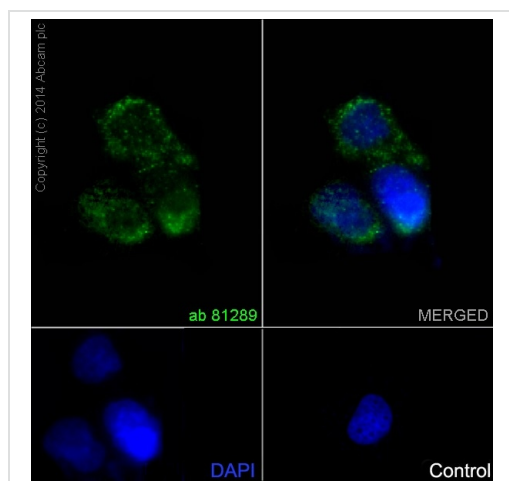
**Lanes 1 - 2:** Merged signal (red and green). Green - Anti-CD59 antibody [EPR6425(2)] (**ab133707**) observed at 14 kDa. Red - loading control, **ab9484**, observed at 37 kDa.

**ab133707** was shown to specifically react with CD59 in wild-type HAP1 cells as signal was lost in CD59 knockout cells. Wild-type and CD59 knockout samples were subjected to SDS-PAGE. Ab133707 and **ab9484** (Mouse anti-GAPDH loading control) were incubated overnight at 4°C at 1/1000 dilution and 1/20000 dilution respectively. Blots were developed with Goat anti-Rabbit IgG H&L (IRDye® 800CW) preabsorbed **ab216773** and Goat anti-Mouse IgG H&L (IRDye® 680RD) preabsorbed **ab216776** secondary antibodies at 1/20000 dilution for 1 hour at room temperature before imaging.



Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti-CD34 antibody [EP373Y]

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) analysis of human angiosarcoma labeling CD34 with unpurified **ab81289** at 1/100-1/250.



Immunocytochemistry/ Immunofluorescence - Anti-CD34 antibody [EP373Y]

Immunocytochemistry/Immunofluorescence analysis of HUVEC (Human umbilical vein endothelial cell line) cells labelling CD34 with purified **ab81289** at 1/100. Cells were fixed with 4% paraformaldehyde and permeabilized with 0.1% Triton X-100. An Alexa Fluor® 488-conjugated goat anti-rabbit IgG (**ab150077**) (1/500) was used as the secondary antibody. DAPI (blue) was used as the nuclear counter stain.

Control: primary antibody (1/100) and secondary antibody, **ab150120**, an Alexa Fluor® 488-conjugated goat anti-mouse IgG (1/500).

Why choose a recombinant antibody?

**Research with confidence**  
Consistent and reproducible results

**Long-term and scalable supply**  
Recombinant technology

**Success from the first experiment**  
Confirmed specificity

**Ethical standards compliant**  
Animal-free production

Hematopoietic Stem Cell Marker (CD34, CD59, CD90 / Thy1, CD38, c-Kit) Antibody Panel - Human (ab254022)

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Promise 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