

# PRODUCT INFORMATION



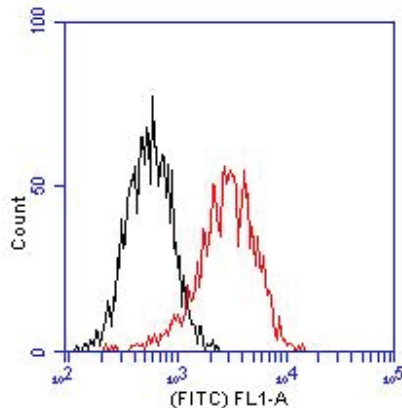
## CD34 Monoclonal FITC Antibody (Clone ICO-115)

Item No. 10208

### Overview and Properties

**Contents:** This vial contains 500 µg of ammonium sulfate-purified, fluorescein labeled anti-CD34 IgG1  
**Synonym:** Hematopoietic Progenitor Cell Antigen CD34  
**Cross Reactivity:** (+) CD34  
**Species Reactivity:** (+) Human  
**Uniprot No.:** P28906  
**Form:** Liquid  
**Storage:** 4°C (as supplied)  
**Stability:** ≥6 months  
**Storage Buffer:** PBS, pH 7.2  
**Clone:** ICO-115  
**Host:** Mouse  
**Applications:** Flow cytometry (FC) and Immunofluorescence (IF); the recommended starting dilution is 20-50 µg/ml for FC and IF. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

### Image



Flow cytometry of human erythroleukemia cells.  
Peak 1: IgG1-FITC isotype control  
Peak 2: CD34 FITC (50 µg/ml)

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/08/2021

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

CD34 is a transmembrane phosphoglycoprotein and sialomucin protein that is commonly used as a marker for hematopoietic progenitor cells.<sup>1,2</sup> It is composed of an N-terminal signaling peptide, a sialylated and O-glycosylated extracellular mucin domain, a cysteine-containing globular domain, a juxtamembrane stalk region, and an intracellular C-terminal tail.<sup>3</sup> CD34 is expressed primarily in hematopoietic progenitor cells but is also expressed in mesenchymal stromal, muscle satellite, interstitial, epithelial and vascular endothelial progenitor, and adipose mesenchymal stem/stromal cells.<sup>1,3</sup> It has both pro- and anti-adhesion activities, providing an anchor for L-selectin-mediated attachment of lymphocytes in vascular endothelial cells and blocking adhesion of bone marrow-derived mast cells *in vitro*.<sup>1,4</sup> CD34 has commonly been used as a marker for the selection and enrichment of hematopoietic stem cells for bone marrow transplants.<sup>1</sup> Cayman's CD34 Monoclonal FITC Antibody (Clone ICO-115) can be used for flow cytometry (FC) and immunofluorescence (IF) applications.

## References

---

1. Sidney, L.E., Branch, M.J., Dunphy, S.E., *et al.* Concise review: Evidence for CD34 as a common marker for diverse progenitors. *Stem Cells* **32(6)**, 1380-1389 (2014).
2. Hughes, M.R., Hernaez, D.C., Cait, J., *et al.* A sticky wicket: Defining molecular functions for CD34 in hematopoietic cells. *Exp. Hematol.* **86**, 1-14 (2020).
3. Scherberich, A., Di Maggio, N., and McNagny, K.M. A familiar stranger: CD34 expression and putative functions in SVF cells of adipose tissue. *World J. Stem Cells* **5(1)**, 1-8 (2013).
4. Drew, E., Merzaban, J.S., Seo, W., *et al.* CD34 and CD43 inhibit mast cell adhesion and are required for optimal mast cell reconstitution. *Immunity* **22(1)**, 43-57 (2005).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM