

# PRODUCT INFORMATION



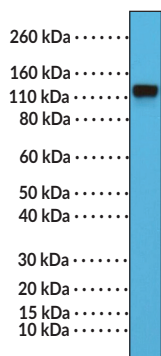
## Integrin $\beta 1$ /CD29 Rabbit Monoclonal Antibody (Clone RM285)

Item No. 32237

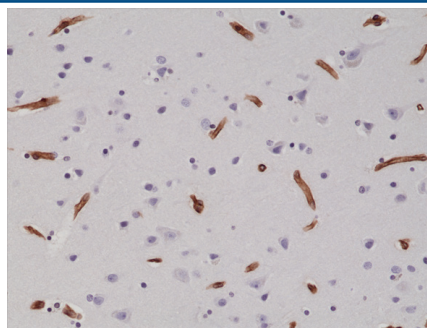
### Overview and Properties

<b>Contents:</b>	This vial contains 100 $\mu$ l of protein A-affinity purified monoclonal antibody.
<b>Synonyms:</b>	Cluster of Differentiation 29, GPIIa, ITGB1, Fibronectin Receptor Subunit $\beta$ , Platelet Glycoprotein IIa, VLA-4 Subunit $\beta$
<b>Immunogen:</b>	Peptide corresponding to human integrin $\beta 1$
<b>Cross Reactivity:</b>	(+) Integrin $\beta 1$
<b>Species Reactivity:</b>	(+) Human
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	$\geq 1$ year
<b>Storage Buffer:</b>	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Clone:</b>	RM285
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 for IHC and 1:2,000-1:4,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

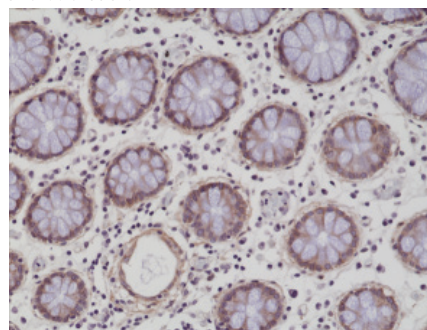
### Images



WB of HeLa cell lysates using Integrin  $\beta 1$ /CD29 Rabbit Monoclonal Antibody (Clone RM285) at a dilution of 1:4,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human brain tissue using Integrin  $\beta 1$ /CD29 Rabbit Monoclonal Antibody (Clone RM285) at a 1:200 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human colon tissue using Integrin  $\beta 1$ /CD29 Rabbit Monoclonal Antibody (Clone RM285) at a 1:200 dilution.

**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, 01/30/2024

**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

---

Integrins are cell adhesion receptors that mediate the attachment of cells to the extracellular matrix (ECM) and facilitate cell-cell interactions.<sup>1</sup> The integrin family contains 18  $\alpha$  subunits and eight  $\beta$  subunits that form 24 different heterodimers grouped into familial subgroups based on subunit composition and/or their ligand binding properties. Integrin  $\beta 1$ , also known as CD29, is an integrin subunit that heterodimerizes with  $\alpha$  integrin subunits, forming 12 different  $\alpha$ - $\beta$  integrin heterodimers that function as receptors for a variety of ECM proteins, including collagens, fibronectins, and laminins, and mediate outside-in and inside-out signaling cascades that regulate numerous cellular processes, including proliferation, adhesion, survival, migration, and organogenesis.<sup>2</sup> Integrin  $\beta 1$  is expressed in most tissues, with high levels found in epidermal stem cells.<sup>1,3</sup> Increased tumor levels of integrin  $\beta 1$  have been found in patients with lung or breast cancer and are associated with decreased overall survival.<sup>4</sup> Cayman's Integrin  $\beta 1$ /CD29 Rabbit Monoclonal Antibody (Clone RM285) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

## References

---

1. Barczyk, M., Carracedo, S., and Gullberg, D. Integrins. *Cell Tissue Res.* **339(1)**, 269-280 (2010).
2. Howe, G.A. and Addison, C.L.  $\beta 1$  integrin: An emerging player in the modulation of tumorigenesis and response to therapy. *Cell Adh. Migr.* **6(2)**, 71-77 (2012).
3. Brakebusch, C., Hirsch, E., Potocnik, A., *et al.* Genetic analysis of  $\beta 1$  integrin function: Confirmed, new and revised roles for a crucial family of cell adhesion molecules. *J. Cell Sci.* **110(Pt 23)**, 2895-2904 (1997).
4. Sun, Q., Zhou, C., Ma, R., *et al.* Prognostic value of increased integrin-beta 1 expression in solid cancers: A meta-analysis. *Onco Targets Ther.* **11**, 1787-1799 (2018).

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