



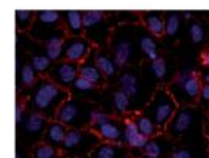
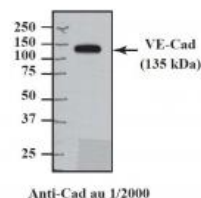
## VE-Cadherin Anti-Cad 3 Polyclonal Antibody

Cat No: A03144 - 100 µg

### General Data

<b>Shipping:</b>	Dry ice
<b>Stability:</b>	Store at -20degrees;shelf life 2 years maximum after production
<b>Formulation:</b>	liquid
<b>Storage Buffer:</b>	0.1M Tris Glycine pH 7.8, 0.02% Na azide
<b>Host:</b>	Rabbit
<b>Antigen:</b>	Fragment of mature VE-Cadherin

**Application(s):** WB, IF



Anti-Cad au 1/2000

Anti-Cad3 au 1/200

### Product Overview

Cadherins are calcium-dependent cell adhesion proteins. They preferentially interact with themselves in a homophilic manner in connecting cells; cadherins may thus contribute to the sorting of heterogeneous cell types. VE-cadherin may play an important role in endothelial cell biology through control of the cohesion and organization of the intercellular junctions. It associates with alpha-catenin forming a link to the cytoskeleton. It acts in concert with KRIT1 to establish and maintain correct endothelial cell polarity and vascular lumen.

### Scientific Literature

Hermant B, Bibert S, Concord E, Dublet B, Weidenhaupt M, Vernet T, Gulino-Debrac D. Identification of proteases involved in the proteolysis of vascular endothelium cadherin during neutrophil transmigration. *J Biol Chem.* 2003 Apr 18;278(16):14002-12.

Gulin

FP/22/20

For research laboratory use only – Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

#### Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac  
10 bis avenue Ampère  
78180 Montigny le Bretonneux - France  
Tel.: +33 (0)139 306 036

<https://www.bertin-bioreagent.com/pa206/contact-us>  
<https://www.bertin-bioreagent.com>