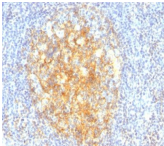


CD106 / VCAM-1 Monoclonal Mouse Antibody (VCAM1/843)



Product Description

CD106 (also known as vascular cell adhesion molecule-1, or VCAM-1) is a member of the Ig superfamily of adhesion molecules. It is expressed at high levels on cytokine stimulated vascular endothelial cells, and at minimal levels on un-stimulated endothelial cells. It is also present on follicular and inter-follicular dendritic cells of lymph nodes, myoblasts, and some macrophages. CD106 serves as a ligand for leukocyte beta-1 integrin VLA-4 and mediates cell adhesion of leukocytes to activated endothelium. It plays a role in various immunological and inflammatory responses. Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 0843, Anti-CD106 (VCAM1/843)**

| Product attributes | |
|---------------------------------------|--|
| Antibody number | #0843 |
| Antibody reactivity (target) | CD106, VCAM-1 |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal |
| Clone | VCAM1/843 |
| Isotype | IgG1, kappa |
| Molecular weight | ~110 kDa |
| Synonyms | CD106; VCAM-1; Vascular Cell Adhesion Molecule 1; VCAM1; INCAM-100 |
| Human gene symbol | VCAM1 |
| Entrez gene ID | 7412 |
| SwissProt | P19320 |
| Unigene | 109225 |
| Immunogen | Recombinant human VCAM1 protein |
| Verified antibody applications | IHC (FFPE) (verified) |
| Antibody target cellular localization | Plasma membrane |
| Species reactivity | Human |
| Antibody application notes | Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunofluorescence: 1-2 ug/mL. Immunohistology formalin-fixed 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Optimal dilution for a specific application should be determined by user |
| Positive control | Human placenta or tonsil |
| Shipping condition | Room temperature |
| Storage Conditions | Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended |
| Regulatory status | For research use only (RUO) |
| Antibody/conjugate formulation | Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide. HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA. Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide. Purified, BSA-free: 1 mg/mL in PBS without azide |
| Antibody research areas | Immunology |
| Cell/tissue expression | Endothelial cells |

| Antibody # prefix | Conjugation | Ex/Em (nm) | Laser line | Detection channel |
|-------------------|--------------------|------------|------------|--------------------------|
| BNC04 | CF®405S | 404/431 | 405 | DAPI (microscopy), AF405 |
| BNC88 | CF®488A | 490/515 | 488 | GFP, FITC |
| BNC68 | CF®568 | 562/583 | 532, 561 | RFP, TRITC |
| BNC94 | CF®594 | 593/614 | 561 | Texas Red® |
| BNC40 | CF®640R | 642/662 | 633-640 | Cy®5 |
| BNC47 | CF®647 | 650/665 | 633-640 | Cy®5 |
| BNCB | Biotin | N/A | N/A | N/A |
| BNUB | Purified | N/A | N/A | N/A |
| BNUM | Purified, BSA-free | N/A | N/A | N/A |

Dye Features
[CF®405S Features](#)
[CF®488A Features](#)
[CF®568 Features](#)
[CF®594 Features](#)
[CF®640R Features](#)
[CF®647 Features](#)

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, LI-COR Bioscience.