

# PRODUCT INFORMATION

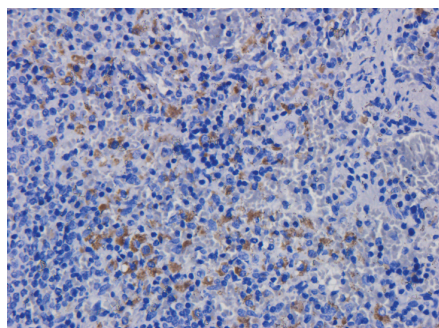


## ITGAL/CD11a Chimeric Monoclonal Antibody (Clone hu1124 (Efalizumab)) Item No. 37155

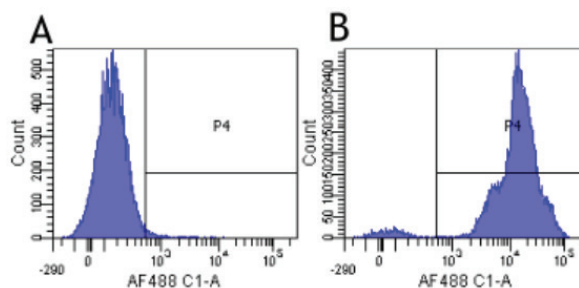
### Overview and Properties

<b>Contents:</b>	This vial contains 200 µg of protein A-affinity purified monoclonal antibody.
<b>Synonyms:</b>	CD11 Antigen-like Family Member A, Integrin subunit αL, ITGAL, Leukocyte Adhesion Glycoprotein LFA-1 α chain, LFA-1A
<b>Immunogen:</b>	Recombinant human CD11a
<b>Species Reactivity:</b>	(+) Human; other species not tested
<b>Uniprot No.:</b>	P20701
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS with 0.02% ProClin™ 300
<b>Concentration:</b>	1 mg/ml
<b>Clone:</b>	hu1124 (Efalizumab)
<b>Host:</b>	Chimeric Monoclonal Antibody
<b>Isotype:</b>	IgG1k
<b>Applications:</b>	Flow Cytometry (FC); block, working concentration/dilution should be determined empirically.

### Images



Immunohistochemical staining of formalin-fixed and paraffin-embedded rat spleen tissue using ITGAL/CD11a Chimeric Monoclonal Antibody (Clone hu1124 (Efalizumab)) at a concentration of 3 µg/ml.



Flow cytometry using the ITGAL/CD11a Chimeric Monoclonal Antibody (Clone hu1124 (Efalizumab)) (Item No. 37155). Human lymphocytes were stained with an isotype control (panel A) or the rabbit-chimeric version of Efalizumab (panel B) at a concentration of 1 µg/ml for 30 minutes at RT. After washing, bound antibody was detected using a AF488 conjugated donkey anti-rabbit antibody and cells analysed on a FACSCanto flow-cytometer.

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, 02/13/2024

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

---

CD11a, also known as integrin  $\alpha$ L (ITGAL), is a protein subunit of the heterodimeric leukocyte integrin  $\alpha$ L $\beta$ 2.<sup>1</sup> It is composed of an extracellular domain that contains an I-domain and a metal-binding domain, which are important for binding, a transmembrane region, and a short cytoplasmic tail.<sup>2</sup> CD11a is expressed ubiquitously on the surface of leukocytes as a component of  $\alpha$ L $\beta$ 2, which binds to intracellular adhesion molecule 1-5 (ICAM-1-5), junctional adhesion molecule A (JAM-A), and endothelial cell-specific molecule 1 (ESM-1) to facilitate T cell adhesion and migration during extravasation into the lymph nodes and tissue.<sup>3</sup> SNPs in *ITGAL* are associated with a reduced susceptibility to Behçet's disease, a systemic inflammatory disease.<sup>4</sup> Cayman's ITGAL/CD11a Chimeric Monoclonal Antibody (Clone hu1124 (Efalizumab)) was recombinantly generated based on the sequence of the original humanized anti-CD11a antibody hu1124.<sup>5</sup> It can be used for flow cytometry (FC).

## References

---

1. Bednarczyk, M., Stege, H., Grabbe, S., *et al.*  $\beta$ 2 Integrins—multi-functional leukocyte receptors in health and disease. *Int. J. Mol. Sci.* **21**(4), 1402 (2020).
2. Knight, J.M., Lee, S.-H., Roberts, L., *et al.* CD11a polymorphisms regulate TH2 cell homing and TH2-related disease. *J. Allergy Clin. Immunol.* **133**(1), 189-197 (2014).
3. Walling, B.L. and Kim, M. LFA-1 in T cell migration and differentiation. *Front. Immunol.* **9**, 952 (2018).
4. Park, S.R., Park, K.S., Park, Y.J., *et al.* CD11a, CD11c, and CD18 gene polymorphisms and susceptibility to Behçet's disease in Koreans. *Tissue Antigens* **84**(4), 398-404 (2014).
5. Bauer, R.J., Dedrick, R.L., White, M.L., *et al.* Population pharmacokinetics and pharmacodynamics of the anti-CD11a antibody hu1124 in human subjects with psoriasis. *J. Pharmacokinet. Biopharm.* **27**(4), 397-420 (1999).

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