

MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT α V β 1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C3001-1

Lot Number: C3001-1-45B-071717

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Cellbanker 2

Host cell: CHO B2

Transfection: Expression vector containing full-length human ITGAV cDNA (GenBank accession number NM_002210.4) and full-length human ITGB1 (GenBank accession number NM_002211.3)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS, 10 μ g/mL puromycin, 250 μ g/mL hygromycin

Stability: Stable for a minimum of 2 months in continuous culture

Background: α V β 1 integrins are transmembrane heterodimeric receptors that contribute to diverse biological functions and play critical roles in many human diseases. α V β 1 is highly expressed on activated fibroblasts, directly binds to the latency-associated peptide of transforming growth factor- β 1 (TGF β 1) and mediates TGF β 1 activation.

Application:

Figure 1

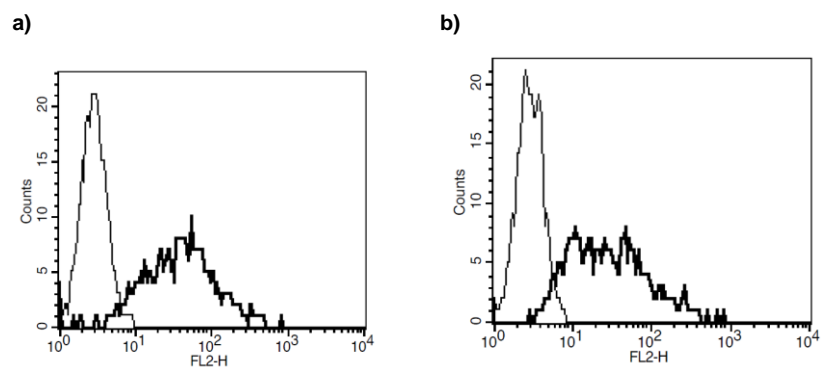


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using using CD51 antibody (a) and CD29 antibody (b). Thin line: parental cells; thick line: receptor-expressing cells.

References:

Nilgun I. Reed, et al. (2003) The avb1 integrin plays a critical in vivo role in tissue fibrosis. *Science Translational Medicine*. 20 May 2015 Vol 7 Issue 288 288ra79

FOR RESEARCH USE ONLY.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.