



# Mouse (monoclonal) Anti-Human CD44 variant (v6) Biotin Conjugate

## PRODUCT ANALYSIS SHEET

<b>Catalog Number:</b>	AHS4469
<b>Lot Number:</b>	See product label
<b>Clone Number:</b>	VFF-7
<b>Quantity/Volume:</b>	100 tests/1 mL
<b>Form of Antibody:</b>	Biotin conjugated purified immunoglobulin in 50 mM Tris, 100 mM NaCl, pH 7.4, with 1% BSA and 0.02% sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
<b>Purity:</b>	Greater than 95% as determined by SDS gel electrophoresis prior to the addition of carrier protein and preservative.
<b>Isotype:</b>	IgG1
<b>Specificity:</b>	<p>CD44 is the receptor for hyaluronan. The interaction between CD44 and hyaluronan may mediate binding of cells to basement membranes. Several CD44 isoforms or variants arise from alternative splicing of mRNA transcribed from a single gene located on chromosome 11. The alternatively spliced mRNA's are translated into a family of protein isoforms which contain inserts in their cytoplasmic domains of varying lengths.</p> <p>Most cell types express the standard CD44, while the expression of the isoforms resulting from the alternative splicing is of limited distribution. CD44 isoforms resulting from alternative splicing have been detected in many human tumors; therefore, the identification of CD44 isoforms has utility in cancer studies.</p> <p>The epitope recognized by this antibody is a variant portion of CD44 encoded by exon v6. While the standard CD44 is observed to have ubiquitous expression, CD44 v6 expression appears to be limited to activated lymphocytes and macrophages.</p>
<b>Applications:</b>	<p>This antibody is suitable for use in flow cytometry. Use 10 µL to label 10<sup>6</sup> cells. This antibody is also suitable for use in immunohistochemistry with cryostat sections, paraffin-embedded sections (after microwave treatment), and cell smears, and Western blotting. The optimal antibody concentration should be determined for each specific application.</p>

**This product is for research use only. Not for use in diagnostic procedures.**

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI AHS4469

(Rev 11/08) DCC-08-1089

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

**Storage:** Store at 2-8°C.

**Expiration Date:** Expires one year from date of receipt when stored as instructed.

**References:**

Arch, R., K. Wirth, M. Hofmann, H. Ponta, S. Matzku, P. Herrlich, and M. Zöller (1992) Participation in normal immune responses of a metastasis-inducing splice variant of CD44. *Science* 257:682-685.

Günthert, U., M. Hofmann, W. Rudy, S. Reber, M. Zöller, I. Haußmann, S. Matzku, A. Wenzel, H. Ponta, and P. Herrlich (1991) A new variant of glycoprotein CD44 confers metastatic potential to rat carcinoma cells. *Cell* 65:13-24.

Heider K.-H., M. Hofmann, E. Horst, F. van den Berg, H. Ponta, P. Herrlich, and S. Pals (1993) A human homologue of the rat metastasis-associated variant of CD44 is expressed in colorectal carcinomas and adenomatous polyps. *J. Cell. Biol.* 120:227-233.

Herrlich, P., W. Rudy, M. Hofmann, R. Arch, M. Zöller, V. Zawadzki, C. Tölg, A. Hekele, S. Pals, K.-H. Heider, J. Sleeman, and H. Ponta (1993) CD44 and splice variants in normal differentiation and tumor progression. In "Cell adhesion molecules". Eds.: D.M. Livingston et al., Edigraf, Rome.

Hofmann, M., W. Rudy, M. Zöller, C. Tölg, H. Ponta, P. Herrlich, and U. Günthert (1991) CD44 splice variants confer metastatic behavior in rats: Homologous sequences are expressed in human tumor cell lines. *Cancer Res.* 51:5292-5297.

Koopman, G., K.-H. Heider, E. Horst, G.R. Adolf, F.v.d. Berg, H. Ponta, P. Herrlich, and S.T. Pals (1993) Activated human lymphocytes and aggressive non-Hodgkin's lymphomas express a homologue of the rat metastasis-associated variant of CD44. *J. Exp. Med.* 177:897-904.

Matsumura, Y., and D. Tarin (1992) Significance of CD44 gene products for cancer diagnosis and disease evaluation. *Lancet* 340:1053-1058.

Stamenkovic, I., A. Aruffo, M. Amiot, and B. Seed (1991) The hematopoietic and epithelial forms of CD44 are distinct polypeptides with different adhesion potentials for hyaluronate-bearing cells. *EMBO J.* 10:343-348.

Underhill, C. (1992) CD44: The hyaluronan receptor. *J. Cell Science* 103:293-298.

**This product is for research use only. Not for use in diagnostic procedures.**

[www.invitrogen.com](http://www.invitrogen.com)

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: [techsupport@invitrogen.com](mailto:techsupport@invitrogen.com)

PI AHS4469

(Rev 11/08) DCC-08-1089

**Important Licensing Information** - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, [www.invitrogen.com](http://www.invitrogen.com)). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.