



Mouse (monoclonal) Anti-Monkey CD3 FITC Conjugate

PRODUCT ANALYSIS SHEET

Catalog Number:	APS0308
Lot Number:	See product label
Expiration Date:	See product label
Clone:	FN-18
Size:	100 tests, lyophilized
Formulation:	Fluorescein isothiocyanate conjugated purified immunoglobulin, lyophilized.
Buffer:	Prior to lyophilization: 0.5 mL PBS + 1% BSA + 125 mM trehalose
Purification:	Purified from tissue culture supernatant by ion exchange chromatography. Sterile filtered (0.2 µm).
Myeloma/Fusion Partners:	Mouse spleen cells were fused with the SP2/0 Ag14 mouse myeloma cell line.
Immunogen:	Rhesus monkey peripheral blood leukocytes.
Isotype:	Mouse IgG ₁
Specificity:	The antigen recognized by this monoclonal antibody is the rhesus monkey RhT3 antigen (homologue of the human CD3 antigen) with a molecular mass of 20 to 25 kDa. This monoclonal antibody reacts with CD3 from <i>Macaca mulatta</i> (rhesus monkey), <i>Macaca arctoides</i> (stumptail monkey), <i>Macaca fascicularis</i> (cynomolgus monkey), <i>Cercocebus atys</i> (sooty mangabey), <i>Cercopithecus aethiops</i> (African green monkey), and <i>Papio anubis</i> (baboon). This antibody does not react with <i>Callithrix jacobus</i> (marmoset), nor does this antibody react with any of the Hominoidea tested (human, chimpanzee, gorilla, and orangutan).
Reconstitution:	Reconstitute with 0.5 mL sterile distilled water. Further dilute to 5 mL with phosphate buffered saline containing 1% BSA for the working dilution. The addition of 0.02% sodium azide is recommended as a preservative.
Application:	This antibody is suitable for use in flow cytometry. Use 20 to 50 µL of the working dilution to label 10 ⁶ cells (PBMCs).
Storage:	Lyophilized product is stable for at least one year at 2-8°C. After reconstitution, the contents can be safely stored at 2-8°C for one month or for one year at -20°C. Add 0.02% sodium azide to prevent bacterial growth.

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

www.invitrogen.com

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

PIAPS0308

(Rev 03/10) DCC-10-0556

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). 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.

Mouse (monoclonal) Anti-Monkey CD3 FITC Conjugate

References:

- Amara, R.R., Villinger, F., et al. (2001) Control of a mucosal challenge and prevention of AIDS by a multiprotein DNA/MVA vaccine. *Science* 292(5514):69-74.
- Ansari, A.A., et al. (2002) Administration of recombinant rhesus interleukin-12 during acute simian immunodeficiency virus (SIV) infection leads to decreased viral loads associated with prolonged survival in SIVmac251-infected rhesus macaques. *J. Virology* 76(4):1731-1743.
- Barouch, D.H., et al. (2000) Control of viremia and prevention of clinical AIDS in rhesus monkey by cytokine-augmented DNA vaccination. *Science* 290:486-492.
- Brice, G.T., et al. (2000) A novel role for tumor necrosis factor- α in regulating susceptibility of activated CD4+ T cells from human and nonhuman primates for distinct coreceptor using lentiviruses. *Journal of Acquired Immune Deficiency Syndromes* 24(1):10-22 (cites the use of monoclonal antibody FN-18 coated onto tosyl activated coated beads for cell stimulation).
- Chakrabarti, L.A., et al. (2000) Age-dependent changes in T-cell homeostasis and SIV load in sooty mangabeys. *J. Med. Primatol.* 29:158-165.
- Chakrabarti, L.A., et al. (2000) Normal T-cell turnover in sooty mangabeys harboring active Simian Immunodeficiency Virus infection. *J. Virol.* February 2000:11209-11223.
- Croix, D.A., et al. (2000) Effect of myobacterial infection on virus loads and disease progression in Simian Immunodeficiency Virus-infected rhesus monkey. *AIDS Research and Human Retroviruses* 16(17):1895-1908.
- Earl, P.L., et al. (2002) Comparison of vaccine strategies using recombinant env-gag-pol MVA with or without an oligomeric env protein boost in the SHIV rhesus macaque model. *Virology* 294(2):270-281.
- Giavedoni, L.D., et al. (2001) Expression of the interleukin-18 gene from rhesus macaque by the Simian Immunodeficiency Virus does not result in increased viral replication. *J. Interferon Cytokine Res.* 21:173-180.
- Harouse, J.M., et al. (2001) Mucosal transmission and induction of simian AIDS by CCR5-specific simian/human immunodeficiency virus SHIVSF162P3. *J. Virol.* 75(4):1990-1995.
- Keller, E.T., et al. (2001) The impact of chronic estrogen deprivation on immunologic parameters in the ovariectomized rhesus monkey (*Macaca mulatta*) model of menopause. *J. Reprod. Immunol.* 50(1):41-55.
- Kou, Z.-C., et al. (1998) In vivo effects of a bacterial superantigen on macaque TCR repertoires. *J. Immunol.* 160:5170-5180.
- Lafont, B.A.P., et al. (2000) One-round determination of seven leukocyte subsets in rhesus macaque blood by flow cytometry. *Cytometry* 41:193-202.
- Locher, C.P., et al. (2003) Expression patterns of phenotypic markers on lymphocytes from human immunodeficiency virus type 2-infected baboons. *AIDS Research and Human Retroviruses* 19(1):31-40.
- Ma, S., et al. (1997) Genetic construction and characterization of an anti-monkey CD3 single-chain immunotoxin with a truncated diphtheria toxin. *Bioconjug. Chem.* 8(5):695-701.
- Mohri, H., et al. (1998) Rapid turnover of T lymphocytes in SIV-infected rhesus macaques. *Science* 279(5354):1223-1227.
- Neville, D.M. Jr, et al. (1996) A new reagent for the induction of T-cell depletion, anti-CD3-CRM9. *J. Immunother. Emphasis Tumor Immunol.* 19(2):85-92.
- Nooij, F.J., et al. (1986) Differentiation antigens on rhesus monkey lymphocytes. II. Characterization of RhT3, a CD-3-like antigen on T cells. *Eur. J. Immunol.* 16(8):981-984.
- Nooij, F.J., et al. (1986) Polymorphism for RhT3, a CD3-like cell surface antigen, expressed on rhesus monkey T lymphocytes. *Immunology* 59(4):611-620.

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

www.invitrogen.com

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

PIAPS0308

(Rev 03/10) DCC-10-0556

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). 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.

Mouse (monoclonal) Anti-Monkey CD3 FITC Conjugate

Schramm, B., et al. (2000) Cytopathicity of human immunodeficiency virus type 2 (HIV-2) in human lymphoid tissue is coreceptor dependent and comparable to that of HIV-1. *J. Virol.* Oct.: 9594-9600.








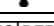

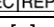

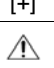
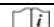
Teranishi, K., et al. (2002) Depletion of anti-gal antibodies in baboons by intravenous therapy with bovine serum albumin conjugated to gal oligosaccharides. *Transplantation* 73(1):129-139.

Thomas, J.M., et al. (1997) Preclinical studies of allograft tolerance in rhesus monkeys: a novel anti-CD3-immunotoxin given peritransplant with donor bone marrow induces operational tolerance to kidney allografts. *Transplantation* 64(1):124-135.

Wilson, L.A., et al. (2000) Identification of SIV env-specific CTL in the jejunal mucosa in vaginally exposed, seronegative rhesus macaques (*Macaca mulatta*). *J. Med. Primatol.* 29:173-181.

Xu, H., et al. (2003) Studies investigating pretransplant donor-specific blood transfusion, rapamycin, and the CD154-specific antibody IDEC-131 in a nonhuman primate model of skin allotransplantation. *J. Immunol.* 170(5):2776-2782.

Explanation of symbols

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		<i>In vitro</i> diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

For Research Use Only. CAUTION: Not for human or animal therapeutic or diagnostic use.

www.invitrogen.com

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

PIAPS0308

(Rev 03/10) DCC-10-0556

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). 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.