

ImmunoTox three-color combinations: iTox3

Immunotoxicology Reagents Now Available
for Non-Human Primate Lymphocyte Phenotyping

General Chemistry
Immunodiagnosics
Molecular Diagnostics
Hematology
Centrifugation
Disease Management
Information Systems
Lab Automation
Hemostasis
Flow Cytometry
Primary Care



BR-10907A

iTox3 System

Lymphocyte Immunophenotyping

The Beckman Coulter anti-non human primate (NHP) iTox3 System is a standardized and comprehensive flow cytometry application to address lymphocyte subset immunophenotyping in monkeys.

We have simplified the preparation steps with easy procedures adapted to several different biological samples, with any lymphocyte sub-family level you want to reach. The iTox3 System offers the flexibility to automate the process or to perform your testing manually. Both methods provide the same high level of confidence in results that you count on from Beckman Coulter.

Delivering Standardized Methodology

The Immunotox iTox3 Combinations have been carefully standardized and validated for accurate and reproducible results. Lysing with the VersaLyse™ "Fix-and-Lyse" no wash procedure allows tremendous discrimination with different specimen types.

Smart iTox 3 Immunophenotyping Reagents

Anti-NHP CD3-FITC / CD4-PC7 / CD8-APC and Anti-NHP CD3-FITC / CD20-PC7 / CD16-APC

The anti-NHP iTox3 System features two multicolor reagents. CD3-FITC / CD4-PC7 / CD8-APC targets the T-cell populations as well as the CD4 and CD8 subsets while CD3-FITC / CD20-PC7 / CD16-APC targets the B lymphocytes and the NK cells. The use of non-overlapping fluorochrome conjugated monoclonal antibodies allows the use of PE-conjugated drop-ins and/or a viability dye like the 7-AAD.

- Low-compensation three-color configuration (FITC/PC7/APC) for Non Human Primate (NHP) and for Rat
- Optimized for dual laser flow cytometers
- Rapid and accurate enumeration of T-, B-, and NK-cell populations including CD4⁺ and CD8⁺ T-cell sub-populations
- Combine with PE-conjugated esoteric markers
- Compatible with cell-viability assessment (7-AAD)
- No-wash procedure
- Easy to implement manually
- Easy to automate



	NHP	Rat
First combination: <i>T cells and T cell subsets</i>	CD3-FITC FN18 (Mouse IgG1, anti-Rhesus) CD4-PC7 M-T477* (Mouse IgG2a, anti-Human) CD8-APC B9.11* (Mouse IgG1, anti-Human)	CD3-FITC 1F4 (Mouse IgM, anti-Rat) CD4-PC7 OX-38 (Mouse IgG2a, anti-Rat) CD8-APC OX-8 (Mouse IgG1, anti-Rat)
Second combination: <i>T cells, B cells, and NK cells</i>	CD3-FITC FN18 (Mouse IgG1, anti-Rhesus) CD20-PC7 B9E9* (Mouse IgG2a, anti-Human) CD16-APC 3G8* (Mouse IgG1, anti-Human)	CD3-FITC 1F4 (Mouse IgM, anti-Rat) CD45RA-PC7 OX-33 (Mouse IgG1, anti-Rat) CD161a-APC 10/78 (Mouse IgG1, anti-Rat)
Viability Dye	7-AAD	7-AAD
Additional Drop-Ins	Customer choice of PE-Conjugated marker	Customer choice of PE-Conjugated marker

Straightforward, Optimized Procedure

- No dilution, no centrifugation, no wash
- Minimized whole blood specimen volume: 25 µL

- Dispense:** 25 µL iTox3 combination
- Dispense:** 25 µL whole blood
- Vortex**
- Incubate:** 20 minutes at room temperature, protected from light
- Add:** 1 mL VersaLyse "Fix-and-Lyse" solution
- Vortex**
- Incubate:** 10 minutes at room temperature, protected from light
- Analyze:** Immediately or keep at 2-8°C, protected from light, and analyze within 2 hours.

**Get more results
with less sample**

* mAbs cross-react with cynomolgus, rhesus and baboon monkeys

Walk-Away Sample Prep and Analysis

The integrated Biomek[®] NXP and FC 500 MPL can handle your iTox assay

The powerful dual laser FC 500 flow cytometer easily accommodates the Immunotox Three-color (iTox3) Combinations. For deeper investigation into the immune status, the FC 500 Series offers simplified five-color analysis, implementing the easiest and most advanced method of color compensation, the Advanced Digital Compensation (ADC) method.

The Multi-Platform Loader (MPL) option of the FC 500 Series provides the capability to automate sample loading using microtubes (e.g. Micronic, Ref. M32022) arranged on a 96-well rack. It can also take on a 40-tube rack for standard 12 x 75 mm test tubes, as well as plates.

The Biomek NXP sets a new standard for flexible laboratory solutions. It puts every aspect of liquid handling – including pipetting, diluting, dispensing and combining samples with reagents – into a single, small-footprint that's as powerful and flexible as it is efficient and economical.

*Automate your
Immunotox Assays*



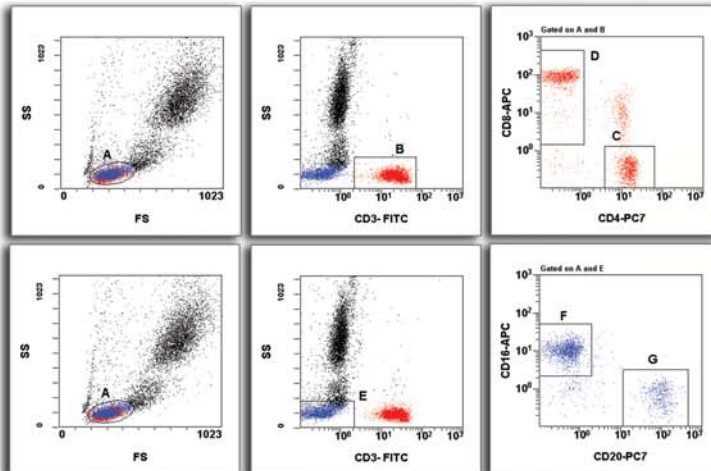
**21 CFR Part 11 Compliance with
both Biomek NXP and FC 500 MPL**

Sensitive and Specific iTox3 Immunophenotyping Reagents

Clear separation between positive and negative events and no background!



Analysis of a Cynomolgus Monkey Whole Blood Sample.

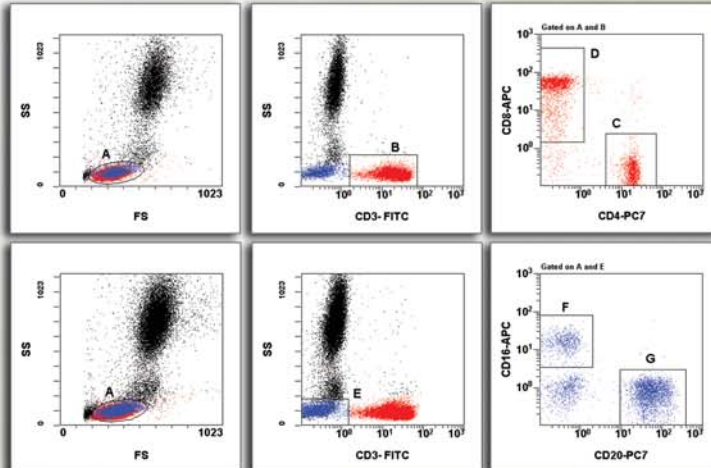


*Get Clear, Accurate
Results, Quickly*

Samples stained with part # A44224 (first row): CD3-FITC, CD4-PC7, and CD8-APC, and with part # A44225 (second row): CD3-FITC, CD20-PC7 and CD16-APC.



Analysis of a Baboon Monkey Whole Blood Sample.



Sample prep with VersaLyse
"Fix-and-Lyse" no wash method.

Ordering Information



	Size	Line	Status	Part No.
Anti-NHP CD3-FITC / CD4-PC7 / CD8-APC	50 tests	IOTest®	RUO	A44224
Anti-NHP CD3-FITC / CD20-PC7 / CD16-APC	50 tests	IOTest	RUO	A44225
Cross-reacting Human Antibodies (single color)				
Anti-Human CD25-RD1 (Clone 1HT44H3)	50 tests	CYTO-STAT®	ASR	6604422
Anti-Human CD69-PE (Clone TP1.55.3)	100 tests	IOTest	ASR	IM1943U
Anti-Human CD86-PE (Clone HA5.2B7)	100 tests	IOTest	ASR	IM2729U



Anti-Rat CD3-FITC / CD4-PC7 / CD8-APC	50 tests	IOTest	RUO	A32909
Anti-Rat CD3-FITC / CD45RA-PC7 / CD161a-APC	50 tests	IOTest	RUO	A32910
Other (single color)				
Anti-Rat CD45-PE (Clone OX-1)	50 tests	IOTest	RUO	A36700

Extend the panel with your choice of conjugated markers from our Custom Design Service.

Viability and Sample Prep Reagents

7-AAD Viability Dye	RUO	IM3422
VersaLyse Lysing Solution	RUO	IM3648
IOTest 3 Fixative Solution	RUO	IM3515

Available System Set-up Reagents

APC (675/633) Set-up Kit	RUO	6607120
Flow-Set™ Fluorospheres	IVD & CE	6607007
Flow-Check™ Fluorospheres	IVD & CE	6605359
Cyto-Trol™ Control Cells Kit	IVD & CE	6604248
Anti-Human CD45-FITC	ASR	IM0782U
Anti-Human CD45-PC7	ASR	IM3548U
Anti-Human CD45-APC	ASR	IM2473U

CAUTION: These reagents are designed for use with certain animal biological samples and are not to be used with human biological samples.
RUO: Research Use Only Reagents • ASR: Analyte Specific Reagents • IVD & CE: In Vitro Diagnostic Products & CE-marked Products

Australia, Gladesville (61) 2 9844-6000 Canada, Mississauga (1) 905 819 1234 China, Beijing (86) 10 6515 6028
Czech Republic, Prague (420) 267 00 85 13 Eastern Europe, Middle East, North Africa, South West Asia: Switzerland, Nyon (41) 22 365 3707
France, Villepinte (33) 1 49 90 90 00 Germany, Krefeld (49) 2151 33 35 Hong Kong (852) 2814 7431
India, Mumbai (91) 22 3080 5101 Italy, Cassina de' Pecchi, Milan (39) 02 953921 Japan, Tokyo (81) 3 5530 8500 Korea, Seoul (82) 2 404 2146
Latin America (1) (305) 380 4709 Mexico, Mexico City (001) 52 55 9183 2800 Netherlands, Mijdrecht (31) 297 230630 Puerto Rico (787) 747 3335
Singapore (65) 6339 3633 South Africa/Sub-Saharan Africa, Johannesburg (27) 11 805 2014/5 Spain, Madrid (34) 91 3836080
Sweden, Bromma (46) 8 564 85 900 Switzerland, Nyon (41) 0800 850 810 Taiwan, Taipei (886) 2 2378 3456 Turkey, Istanbul (90) 216 309 1900
UK, High Wycombe (44) 01494 441181 USA, Fullerton, CA (1) 800 742 2345

