



Monoclonal Mouse Anti-Human HLA-DR Antigen/FITC, Clone AB3
Monoclonal Mouse Anti-Human HLA-DR Antigen/RPE, Clone AB3

Code F7266
Code R7267

Analyte specific reagent. Analytical and performance characteristics are not established.

Summary and explanation HLA-DR antigen is constitutively expressed on antigen-presenting cells, such as B lymphocytes, monocytes and dendritic cells but can also be detected on activated T lymphocytes and activated granulocytes (1, 2). Occasionally, natural killer cells express HLA-DR antigen (3). Anti-Human HLA-DR Antigen, AB3, reacts with the DRw52 determinant of the HLA-DR molecule.

Reagent provided F7266 is a purified monoclonal mouse antibody conjugated with fluorescein isothiocyanate isomer 1 (FITC). R7267 is a purified monoclonal mouse antibody conjugated with R-phycoerythrin (RPE).
The conjugates are provided in liquid form in buffer containing 1% bovine serum albumin (BSA) and 15 mmol/L NaN₃, pH 7.2.
Isotype: IgG2a, kappa. Conjugate concentration mg/L: See label on vial.

Immunogen B cells from a follicular lymphoma.

Precautions

1. Analyte specific reagent. Analytical and performance characteristics are not established.
2. For professional users.
3. This product contains sodium azide (NaN₃), a chemical highly toxic in pure form. At product concentrations, though not classified as hazardous, sodium azide may react with lead and copper plumbing to form highly explosive build-ups of metal azides. Upon disposal, flush with large volumes of water to prevent metal azide build-up in plumbing.
4. Minimize microbial contamination of reagents or increase in non-specific staining may occur.
5. As with any product derived from biological sources, proper handling procedures should be used.
6. Wear appropriate Personal Protective Equipment to avoid contact with eyes and skin.
7. Unused solution should be disposed of according to local, State and Federal regulations.

Storage Store in the dark at 2-8 °C. Do not use after expiration date stamped on vial. If reagents are stored under any conditions other than those specified, the conditions must be verified by the user. There are no obvious signs to indicate instability of this product. If unexpected staining is observed which cannot be explained by variations in laboratory procedures and a problem with the antibody is suspected, contact Dako Technical Services.

References

1. Naverrete CV. The HLA system in blood transfusion. *Baillière's Clin Haematol* 2000;13:511-32.
2. Erber WN, Pinching AJ, Mason DY. Immunocytochemical detection of T and B cell populations in routine blood smears. *Lancet* 1984;i:1042-5.
3. Van Dongen JJM, Adriaansen HJ. Immunobiology of leukaemia. In: Henderson ES, Lister TA, Greaves MF, editors. *Leukemia*. Philadelphia, London, Toronto, Montreal, Sydney, Tokyo: WB Saunders Company; 1996. p. 83-130.

Explanation of symbols

REF	Catalogue number		Keep away from sunlight (consult storage section)		Manufacturer
	Consult instructions for use		Batch code		
	Temperature limitation		Use by		

Manufactured by:
Dako Denmark A/S
Produktionsvej 42
DK-2600 Glostrup

Tel. +45 44 85 95 00
Fax +45 44 85 95 95
www.dako.com

Distributed by:
Dako North America, Inc.
6392 Via Real
Carpinteria, California 93013 USA

Tel. 805 566 6655
Fax 805 566 6688
Technical Support 800 424 0021
Customer Service 800 235 5763