

NOTICE OF DISCONTINUATION

Please note that we will discontinue this item in your region. We will continue to offer the following single-color equivalent replacements, which are labeled Analyte Specific Reagents (ASR) or In Vitro Diagnostics (IVD).

Single-color equivalent replacements:

- F0434 Polyclonal Rabbit Anti-Human Kappa Light Chains/FITC (ASR)
- R0808 Monoclonal Mouse Anti-Human CD19/RPE (IVD)

Learn more about these replacement products by contacting rpsupport@agilent.com or through your local sales representative.

MultiMix™ **Dual-Colour Reagent** **Anti-Human Kappa Light Chains/FITC** **Anti-Human CD19/RPE**

Code FR048

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

Summary and explanation

Most B cells, with the exception of pre-B progenitors, pre-B cells and mature plasma cells, express immunoglobulin on their surface. Each cell expresses only one light chain type. In normal peripheral blood and lymph nodes there is a mixture of kappa-positive and lambda-positive cells, with two-thirds of the cells expressing kappa and one-third expressing lambda. CD19 is the broadest lineage-specific surface marker for B cells and is present on the surface of virtually all B lymphocytes, including early B progenitor cells (1).

Anti-Kappa Light Chains reacts with free kappa chains as well as kappa chains in intact immunoglobulin molecules. Anti-CD19, HD37, was included in the Second, Third, Fourth and Fifth International Workshops and Conferences on Human Leucocyte Differentiation Antigens and studies by a number of laboratories confirmed its reactivity with CD19 (2, 3).

Reagent provided

FR048 comprises the following two, matched, fluorescent antibodies:

Polyclonal Rabbit Anti-Human Kappa Light Chains conjugated with fluorescein isothiocyanate isomer 1 (FITC).

Monoclonal Mouse Anti-Human CD19, Clone HD37, conjugated with R-phycoerythrin (RPE).

The Anti-Kappa Light Chains conjugate has been produced from a F(ab)₂ fragment of an affinity-isolated polyclonal rabbit antibody. The Anti-CD19 conjugate has been produced from a purified monoclonal mouse antibody of isotype IgG1, kappa. FR048 is provided in liquid form in buffer containing 1% bovine serum albumin (BSA) and 15 mmol/L NaN₃, pH 7.2.

Immunogens

Polyclonal Rabbit Anti-Human Kappa Light Chains: Polyclonal immunoglobulin light chains of kappa type isolated from a pool of human sera.

Monoclonal Mouse Anti-Human CD19, Clone HD37: Hairy cell leukemia cells (4).

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. As with any product derived from biological sources, proper handling procedures should be used.
5. Wear appropriate Personal Protective Equipment to avoid contact with eyes and skin.
6. Unused solution should be disposed of according to local, State and Federal regulations.






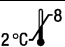

Storage

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 results are observed which cannot be explained by variations in laboratory procedures and a problem with the reagent is suspected, contact Dako Technical Support.

References

1. Pezzutto A, Dörken B, Rabinovitch PS, Ledbetter JA, Moldenhauer G, Clark EA. CD19 monoclonal antibody HD37 inhibits anti-immunoglobulin-induced B cell activation and proliferation. *J Immunol* 1987;138:2793-9.
2. Nadler LM. B cell/leukemia panel workshop: summary and comments. In: Reinherz EL, Haynes BF, Nadler LM, Bernstein ID, editors. *Leukocyte typing II. Proceedings of the 2nd International Workshop on Human Leukocyte Differentiation Antigens*; 1984 Sept 17-20; Boston, USA. New York, Berlin, Heidelberg, Tokyo: Springer-Verlag; 1986. Volume 2. p. 3-43.
3. Mason DY, Ladyman H, Gatter KC. Immunohistochemical analysis of monoclonal anti-B cell antibodies. In: Reinherz EL, Haynes BF, Nadler LM, Bernstein ID, editors. *Leukocyte typing II. Proceedings of the 2nd International Workshop on Human Leukocyte Differentiation Antigens*; 1984 Sept 17-20; Boston, USA. New York, Berlin, Heidelberg, Tokyo: Springer-Verlag; 1986. Volume 2. p. 245-55.
4. Pezzutto A, Dörken B, Feller A, Moldenhauer G, Schwartz R, Wernet P, et al. HD37 monoclonal antibody : a useful reagent for further characterization of "non-T, non-B" lymphoid malignancies. In: Reinherz EL, Haynes BF, Nadler LM, Bernstein ID, editors. *Leukocyte typing II. Proceedings of the 2nd International Workshop on Human Leukocyte Differentiation Antigens*; 1984 Sept 17-20; Boston, USA. New York, Berlin, Heidelberg, Tokyo: Springer-Verlag; 1986. Volume 2. p. 391-402.

Explanation of symbols

 REF	Catalogue number		Keep away from sunlight (consult storage section)		Manufacturer
	Consult instructions for use		LOT		Batch code
	Temperature limitation 2°C - 8°C				Use by

Manufactured by:
Dako Denmark ApS
Produktionsvej 42
DK-2600 Glostrup

Tel. +45 44 85 95 00
Fax +45 44 85 95 95
www.agilent.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

Revision 2020.06