

Datasheet: OBT0009

Description:	RABBIT ANTI HUMAN TdT KIT
Name:	TdT
Other names:	TERMINAL DEOXYNUCLEOTIDYL TRANSFERASE
Format:	Kit
Product Type:	Kits
Isotype:	Polyclonal IgG
Quantity:	20 TESTS

Product Details

Applications

This product has been reported to work in the following applications. This information is derived from testing within our laboratories, peer-reviewed publications or personal communications from the originators. Please refer to references indicated for further information. For general protocol recommendations, please visit www.bio-rad-antibodies.com/protocols.

	Yes	No	Not Determined	Suggested Dilution
Immunofluorescence	■			

Where this product has not been tested for use in a particular technique this does not necessarily exclude its use in such procedures. Suggested working dilutions are given as a guide only. It is recommended that the user titrates the product for use in their own system using appropriate negative/positive controls.

Target Species	Human
Product Form	Kit
Preservative Stabilisers	All components contain 0.1% Sodium Azide (NaN ₃) Components 2, 2C and 3 contain 1% Bovine Serum Albumin Component M contains 80% glycerol
External Database Links	UniProt: P04053 Related reagents Entrez Gene: 1791 DNTT Related reagents
Synonyms	TDT
Product Information	Rabbit anti Human TdT kit is a set of reagents intended to allow the identification of Terminal deoxynucleotidyl transferase positive (TdT ⁺) cells by combined fluorescence and phase contrast microscopy.
Reagents In The Kit	Kit contains ready-to-use reagents sufficient for twenty slide tests: 1 Normal goat serum - 1.0ml

2 Rabbit anti- human -TdT antibody mixture - 0.5ml

2C Negative control antibody - 0.5ml

3 FITC-goat anti- rabbit IgG F(ab')₂ - 1.0ml.

M Mounting medium (80% glycerol in PBS/azide) - 1.0ml

Additional Reagents Required	Cold absolute methanol Phosphate buffered saline (PBS)
---	---

Instructions For Use	Standard bone marrow or peripheral blood smears may be used for immunofluorescence staining. Cytocentrifuge preparations of cell suspensions made up in RPMI-1640 tissue culture medium containing 10% foetal calf serum are also suitable. Frozen sections should be cut as thin as possible. Store unfixed slide preparations in a dessicator at room temperature. 1. Circle two areas of interest (1 cm diameter) with diamond pencil on the slide to be stained. One will be used as a control area and the other is for the test. Fix smears for 30 min. in absolute methanol at 4°C immediately before staining. Air dry. 2. Block by adding a drop of normal goat serum to both circled areas. Let stand for 5 min. Rinse off goat serum with PBS. 3. Apply 25 ul rabbit anti human TdT antibody mixture to the test area and 25 ul of negative control to control area and allow to react for 30 min in a humid chamber. 4. Rinse avoiding cross-contamination of reagents. Now wash for 10 min. in a Coplin jar with several changes of PBS. Agitate occasionally for thorough washing. 5. Wipe the slide, avoiding cells, and shake excess wash solution off circled area. Spread 25 ul of FITC goat anti rabbit IgG F(ab')₂ and incubate in humid chamber for 30 min in the dark. 6. Rinse off and wash for 10 min. in a Coplin jar with several changes of PBS. 7. Apply small drops of glycerol/PBS (mounting medium) to two glass cover slips and invert circled areas on slide over dropped cover slips. 8. Examine by epifluorescence and phase microscopy. Evaluate the percentage of nucleated cells exhibiting nuclear fluorescence.
-----------------------------	---

Storage	Store at +4°C or at -20°C if preferred. Storage in frost-free freezers is not recommended. This product should be stored undiluted. Avoid repeated freezing and thawing as this may denature the antibody. Should this product contain a precipitate we recommend microcentrifugation before use.
----------------	---

Shelf Life	Please see label for expiry date.
-------------------	-----------------------------------

Health And Safety Information	Material Safety Datasheet documentation #10504 #10147 #10143 available at: 1 - Normal goat serum (10504): https://www.bio-rad-antibodies.com/uploads/MSDS/10504.pdf M - Mounting Medium (10147): https://www.bio-rad-antibodies.com/uploads/MSDS/10147.pdf 2, 2C & 3 - Antibody Preparations (10143): https://www.bio-rad-antibodies.com/uploads
--	--

Regulatory

For research purposes only

North & South Tel: +1 800 265 7376

America Fax: +1 919 878 3751

Email: antibody_sales_us@bio-rad.com

Worldwide

Tel: +44 (0)1865 852 700

Fax: +44 (0)1865 852 739

Email: antibody_sales_uk@bio-rad.com

Europe

Tel: +49 (0) 89 8090 95 21

Fax: +49 (0) 89 8090 95 50

Email: antibody_sales_de@bio-rad.com

'M318116:180718'

Printed on 01 Aug 2018