

T-SPOT.® TB Test Results Report

ODL® ID : 991900210

Sample ID : EXAMPLE

00/00/00



Collection Date : 01/11/2023

Received Date : 02/11/2023

Processed On : Received Date

Requesting Laboratory : ZZV10 Validation Order Locaton

Clinician : ZZV10 Dr Single Results

T-SPOT.TB Test Results

Approved By (QC Authoriser) : KLO

Date: 07/11/2023

Assay Result

Positive

A test is Positive if either antigen panel minus the nil control is greater than or equal to 6 spots. A Positive test indicates that Tuberculosis infection is likely.

Nil (Neg) Control Spot Count 0
Panel A Spot Count 15
Panel B Spot Count 15
Positive Control Spot Count >20

Interpretation of Results

- The results should be used and interpreted in the context of the overall clinical picture
 - A full breakdown of interpretation and Quality Control of the test is given at <http://www.oxfordimmunotec.com>
 - A test is positive if either antigen panel (whichever is greater) minus the nil control is greater than or equal to 6 spots. A positive test indicates that Tuberculosis infection is likely
 - A test is negative if antigen panels minus the nil control are less than or equal to 5 spots. A negative test indicates that Tuberculosis infection is unlikely
 - A test is borderline for spot counts around the cutoff (5, 6 or 7 spots) and a retest is recommended. This zone was created as a quality assurance to allow for the variability that can occur with every laboratory test and may be due to potential biological variation. This borderline region represents ODL's uncertainty of measurement for the T-SPOT.TB test
 - A test is indeterminate if the Nil Control spot count is in excess of 10 spots. Another sample should be collected from the individual and tested
 - A test is also considered indeterminate where the Positive Control spot count is <20 spots, unless either Panel A or Panel B is 'Positive' as described in the Results Interpretation and Assay Criteria, in which case the result is valid
- * This result relates only to the whole blood sample received, processed and reported for the T.SPOT.TB assay for the patient ID referenced on this result report

Oxford Diagnostic Laboratories is a trading division of Oxford Immunotec Ltd.
T-SPOT and the Oxford Diagnostic Laboratories logo are registered trademarks of Oxford Immunotec Ltd.
Oxford Immunotec Ltd is trading as Revvity.



Tel: +44 (0) 1235 433164
Fax: +44 (0) 1235 442781
Email : customerexperience@oxfordimmunotec.com
www.oxforddiagnosticlaboratories.eu
Registered in England No. 4516079
VAT No. 201141779

© 2023 Oxford Immunotec Ltd. All rights reserved.

4066