

## Solidscreen II

Microplate for solid phase Antiglobulin Test with TANGO<sup>®</sup> optimo

FOR IN-VITRO DIAGNOSTIC USE

### Package size

[REF] 806521100 [VOL] 10 Micotestplate (12 strips each)

### Intended Use

Solidscreen II is used for the TANGO<sup>®</sup> optimo. The Solidscreen II solid phase antiglobulin test is used as indirect antiglobulin (IAT) test for crossmatch, antibody screening and antibody identification, as well as direct antiglobulin test (DAT) and for the determination of weak D and partial D antigens (DVI and DVII) in donor blood samples.

### Summary

Moreschi first described the use of Anti-Human Globulin in 1908<sup>1</sup>. Coombs rediscovered the test in 1945.<sup>2,3</sup> By injecting rabbits with human IgG, they were able to produce a protein (Anti-IgG) that reacted with incomplete antibodies (IgG). Most "incomplete" antibodies (IgG) fail to agglutinate red blood cells suspended in saline.<sup>4</sup> Most clinically significant antibodies in red blood cell serology are of the IgG class and can only be detected by the use of Anti-IgG. A stable lattice structure is formed and agglutination occurs when Anti-IgG binds to the IgG sensitized red blood cells. The ability to detect alloantibodies or autoantibodies directed against human red blood cells, in human plasma or serum, is a necessary part of routine laboratory testing as well as the detection of weak D and partial D antigens (DVI and DVII) in donors. There are two very important applications for antibody detection:

1. The detection of red blood cell antibodies prior to red blood cell or whole blood transfusion to prevent the possibility of a transfusion reaction with accompanying red cell destruction.
2. To detect the presence of red blood cell antibodies in maternal or newborn serum that may result in Hemolytic Disease of the Newborn.

Routine pretransfusion studies always include tests for alloantibodies or autoantibodies directed against human red blood cells

Routine pretransfusion studies always include tests for the D antigen.

### Principle

Solidscreen II is a solid phase assay for

- a) the detection of red blood cell alloantibodies or autoantibodies in human plasma or serum.
- b) the determination of weak D and partial D antigens (DVI and DVII) of samples which have tested negative with IgM anti-D using Erytype S and the TANGO<sup>®</sup> optimo.

The Solidscreen II well is coated with Protein A. Protein A is a component of the cell wall of *Staphylococcus aureus* and has a very high affinity for the Fc portion of most immunoglobulin classes<sup>5</sup>.

For a) The plasma or serum and Reagent Red Blood Cells are added to the Protein A coated well. Sensitization of the red cell occurs if the corresponding antibody is present for the antigen on the red cell.

For b) Solidscreen II Anti-D Blend Blood Grouping Reagent and test red blood cells are added to the Protein A coated well. Sensitization of the red blood cell occurs if D antigen is present on the red blood cell.

Following incubation, and two wash processes to remove unbound protein, Anti-Human Globulin is added to the well and acts as a link between the antibody coating of neighbouring red blood cells and induces solid phase. Uncoated red blood cells will form a red blood cell button. Following centrifugation, the well is evaluated. A smooth monolayer of red blood cells is indicative of a positive reaction. A compact button of cells in the middle of the well is indicative of a negative reaction.

### Reagent

The Solidscreen II microplate consists of twelve strips containing eight wells per strip. Each well is coated with Protein A. Each Solidscreen II microplate is packaged in a foil container to prevent contamination. Each plate is ready to use.

### Precautions

- For in vitro diagnostic use
- Plates that have been opened and not loaded on the TANGO<sup>®</sup> optimo may be stored, uncovered, in a dry area, not to exceed 24 hours.
- Resuspend Reagent Red Blood Cells prior to use and insert cell mixers before loading on TANGO<sup>®</sup> optimo.
- Do not use beyond the expiration date.
- Do not freeze.
- Do not use beyond seven days on the TANGO<sup>®</sup> optimo
- Do not attempt to reuse unused portions of the strip.
- Let plate come to room temperature before opening the foil packet to limit condensation.
- Store foil packets at 2 to 8°C when not in use.
- Do not use samples collected in gel separator tubes.

### Specimen Collection

#### TANGO<sup>®</sup> optimo

#### For antibody screening and antibody identification (Indirect Antiglobulin Test IAT)

Fresh samples of clotted or EDTA anticoagulated whole blood can be used for antibody screening and antibody identification with the indirect antiglobulin test. Samples collected following standard blood sampling guidelines are acceptable. The specimen should be tested as soon as possible after collection. If testing is delayed, EDTA and clotted specimens should be stored at 2 to 8°C. Use of samples older than seven days should be avoided, since antibody reactivity has been shown to decrease in older samples. Stored samples should be allowed to reach room temperature prior to testing. Blood specimens exhibiting gross hemolysis or contamination should not be used

There must be a distinct separation between the cellular and the plasma layer in the sample tube. Samples can be centrifuged or allowed to settle.

For crossmatch (Indirect Antiglobulin Test)

Fresh samples of EDTA or citrate anticoagulated whole blood samples must be used for the crossmatch. Samples collected following standard blood sampling guidelines are acceptable. The specimen should be tested as soon as possible after collection. If testing is delayed, EDTA specimens should be stored at 2 to 8°C, citrated specimens (donor segments) at 1 to 6°C. Use of EDTA anticoagulated samples older than seven days should be avoided, since antibody reactivity has been shown to decrease in older samples. Donor blood stored in citrate anticoagulant at 1 to 6°C may be tested until the expiration date of the donor unit. These red blood cells to be tested must be prepared prior to testing. Refer to instructions in the TANGO<sup>®</sup> optimo Users Guide<sup>6</sup>. Stored samples should be allowed to reach room temperature prior to testing. Blood specimens exhibiting gross hemolysis or contamination should not be used.

There must be a distinct separation between the cellular and the plasma layer in the sample tube. Samples can be centrifuged or allowed to settle.

#### For Direct Antiglobulin Test (DAT)

Fresh samples of EDTA anticoagulated whole blood samples and cord blood samples must be used for the Direct Antiglobulin Test. Samples collected following standard blood sampling guidelines are acceptable. The specimen should be tested as soon as possible after collection. If testing is delayed, EDTA anticoagulated whole blood samples should be stored at 2 to 8°C. Use of samples older than seven days should be avoided unless there is no other alternative since antibody reactivity has been shown to decrease in older samples. Stored samples should be allowed to reach room temperature prior to testing. Blood specimens exhibiting gross hemolysis or contamination should not be used.

There must be a distinct separation between the cellular and the plasma layer in the sample tube. Samples can be centrifuged or allowed to settle.

#### For weak D and partial D antigen typing (Indirect Antiglobulin Test IAT)

Fresh samples of EDTA or citrate anticoagulated whole blood samples must be used for the weak D test. Samples collected following standard blood sampling guidelines are acceptable. The specimen should be tested as soon as possible after collection. If testing is delayed the EDTA anticoagulated samples should be stored at 2 to 8°C. EDTA anticoagulated whole blood samples may be tested for up to seven days following collection. Donor blood stored in citrate anticoagulant at 1 to 6°C may be tested until the expiration date of the donor unit. The red blood cells to be tested must be prepared prior to testing. Refer to instructions in the TANGO<sup>®</sup> optimo Users Guide.<sup>4</sup> Stored samples should be allowed to reach room temperature prior to testing. Blood specimens exhibiting gross hemolysis or contamination should not be used.

There must be a distinct separation between the cellular and the plasma layer in the sample tube. Samples can be centrifuged or allowed to settle.

### Materials

#### Materials Provided

- Solidscreen II microplates

#### Material required but not provided

- TANGO<sup>®</sup> optimo [REF] 848900010
- Isotonic saline
- MLB 2 (modified LISS Biotest) [REF] 805200100
- Biotestcell<sup>®</sup> Pool [REF] 816065100, Biotestcell<sup>®</sup> 1 & 2 [REF] 816014100, Biotestcell<sup>®</sup> 3 [REF] 816085100, Biotestcell<sup>®</sup>-I 8 [REF] 816020100, Biotestcell<sup>®</sup>-I 11 [REF] 816021100
- Search-Cyte<sup>®</sup> Pool, or Search-Cyte Duo<sup>®</sup>, or Search-Cyte<sup>®</sup> Trio for the TANGO<sup>®</sup> optimo
- Donor or patient red blood cells
- Solidscreen II Anti-D (RH1) Blend [REF] 806530100
- Alsevers Solution [REF] 806510100
- Anti-Human Globulin Anti-IgG Solidscreen II [REF] 806516100
- Solidscreen II Control [REF] 806514100
- Solidscreen II Control B [REF] 806519100
- Solidscreen II Negative Control [REF] 806509100
- PBS pH 7.3 ± 0.2
- Centrifuge
- Cell Mixers

### Test Procedure

#### Indirect Antiglobulin Test (IAT)

#### Crossmatch, antibody screening and antibody identification

1. TANGO<sup>®</sup> optimo dispenses 50µL of patient serum/plasma or control reagents into the Solidscreen II well.
2. TANGO<sup>®</sup> optimo prepares a 1% suspension of Reagent Red Blood Cells with MLB 2. An approx 1% suspension of donor red blood cells is prepared with MLB2 and Alsevers Solution.
3. TANGO<sup>®</sup> optimo dispenses 50µL of the Reagent Red Blood Cells or donor red blood cells prepared in (2.) into the well with patient serum/plasma or control reagents.
4. The mixture is incubated for 20 minutes at 37°C.
5. The mixture is centrifuged following incubation.
6. The supernatant is aspirated and the strip is washed twice. Centrifugation follows each wash process.
7. 100µL of Anti-Human Globulin Anti-IgG Solidscreen II is added to each the well and mixed.
8. Centrifugation by TANGO<sup>®</sup> optimo
9. Reaction is evaluated and interpreted by TANGO<sup>®</sup> optimo.

#### Direct Antiglobulin Test (DAT)

1. TANGO<sup>®</sup> optimo prepares an approx 1% suspension of patient or donor red blood cells with MLB 2 and Alsevers Solution.
2. TANGO<sup>®</sup> optimo dispenses 50µL of the patient or donor red blood cells prepared in (1.) into the Solidscreen II well.
3. Following centrifugation, the supernatant is aspirated and the strip is washed twice. Centrifugation follows each wash process.
4. 100µL of Anti-Human Globulin Anti-IgG Solidscreen II is added to each the well and mixed.
5. Centrifugation by TANGO<sup>®</sup> optimo.
6. Reaction is evaluated and interpreted by TANGO<sup>®</sup> optimo.

**FOR REFERENCE USE ONLY: DO NOT USE in place of package inserts provided with each product.**

