

Blood Grouping Reagents

Anti-A (ABO1)

Seraclone® Murine Monoclonal (A003)

Anti-B (ABO2)

Seraclone® Murine Monoclonal (B005)

Anti-A,B (ABO3)

Seraclone® Murine Monoclonal Blend (BS63/85)

FOR IN-VITRO DIAGNOSTIC USE

For Tube Testing

MEETS FDA POTENCY REQUIREMENTS

U.S. License Number: 1798

Package size

| | | | | |
|-----|-----------|-----|------------|----------------------------|
| REF | 801325100 | VOL | 10 x 10 mL | Seraclone® Anti-A (ABO1) |
| REF | 801350100 | VOL | 10 x 10 mL | Seraclone® Anti-B (ABO2) |
| REF | 801375100 | VOL | 10 x 10 mL | Seraclone® Anti-A,B (ABO3) |

Intended Use

For the determination of the A (ABO1), B (ABO2), A,B (ABO2) antigens of red blood cells using the tube test.

Summary

Between 1900 and 1902, Landsteiner and associates discovered the ABO system of red blood cell antigens. The importance of this discovery is the recognition that antibodies are present when the corresponding antigens are lacking. The ABO system is the only blood group system in which the reciprocal antibodies are consistently and predictably present in most people.¹ Due to this reciprocity, an ABO blood type determination is considered valid if serum typing corresponds with the red blood cell antigen grouping.

| ABO group | Incidence (%) in US population ¹ | |
|-----------|---|--------|
| | Whites | Blacks |
| O | 45 | 49 |
| A | 40 | 27 |
| B | 11 | 20 |
| AB | 4 | 4 |

Biotest Anti-A, Anti-B and Anti-A,B Blood Group Reagents are used to test for the presence or absence of the corresponding antigens. Routine pretransfusion studies always include tests for the ABO antigens.

Phenotype Frequency (%)²

| | Caucasians | Blacks | Asians | Mexican |
|------------------|------------|--------|--------|---------|
| A ₁ | 33 | 19 | 27 | 22 |
| A ₂ | 10 | 8 | Rare | 6 |
| B | 9 | 20 | 25 | 13 |
| O | 44 | 49 | 43 | 55 |
| A ₁ B | 3 | 3 | 5 | 4 |
| A ₂ B | 1 | 1 | Rare | Rare |

Principle of the test

The test principle is hemagglutination. The antibodies in Seraclone® Anti-A (ABO1), Seraclone® Anti-B (ABO2), Seraclone® Anti-A,B (ABO3) bind to the corresponding antigen on red blood cells being tested and cause an antigen-antibody reaction visible as red blood cell agglutination. The four ABO blood groups A, B, A,B and O are defined by the presence or absence of A and B antigens on red blood cells. The absence of both A and B antigens defines blood type O. The antigens A and B react with the corresponding antibody in Seraclone® Anti-A (ABO1), Seraclone® Anti-B (ABO2), and Seraclone® Anti-A,B (ABO3).

Reagent

Seraclone® Anti-A (ABO1), Seraclone® Anti-B (ABO2), and Seraclone® Anti-A,B (ABO3) contain as reactive components monoclonal antibodies of the immunoglobulin class IgM.

They are derived from hybridoma cell lines which are created by fusing mouse antibody producing B lymphocytes with mouse myeloma cells and demonstrate consistent specificity and reproducibility characteristic for monoclonal antibodies. Both antibodies derived from a single clone (sister cells of one hybridoma cell) and a mixture of different antibodies derived from several clones are called monoclonal. Antibodies are diluted in a buffered protein solution containing bovine albumine, ethylenediamine tetraacetate (EDTA), and as colorant Patent Blue (Anti-A) or Tartrazin (Anti-B).

| | |
|----------------------------|------------------------------|
| Seraclone® Anti-A (ABO1) | clone A003 (IgM) |
| Seraclone® Anti-B (ABO2) | clone B005 (IgM) |
| Seraclone® Anti-A,B (ABO3) | clones BS 63/BS 85 (IgM/IgM) |

Preservative: 0.1% sodium azide.

Precautions

- For In-vitro diagnostic use.
- Store at 2 to 8°C.
- Do not use beyond the expiration date.
- Do not use if turbid.
- Handle and dispose of reagents as potentially infectious
- Caution: Do not pipette by mouth. The absence of murine viruses has not been determined.
- Caution: This product contains Natural Rubber Latex Which May Cause Allergic Reactions.
- Warning: Contains sodium azide (NaN₃), which may react with lead or copper plumbing to form explosive azides. If discarded in the sink, flush with large amounts of water to prevent the build-up of explosive metal azides.
- The bovine albumin used for the production of this reagent is purchased from BSE-free US sources, Boval Company L.P. in Cleburne, Tx, USA and Millipore in Kankakee, IL, USA.

Specimen Collection

Fresh samples of clotted, EDTA or citrate anticoagulated whole blood collected following general 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, citrated specimens (donor segments) at 1 to 6°C. Blood specimens exhibiting gross hemolysis or contamination should not be used. Clotted samples or those collected in EDTA may be tested within ten days from collection. Donor blood stored in citrate anticoagulant may be tested until the expiration date of the donor unit.

Materials

Materials provided

- Seraclone® Anti-A (ABO1), Seraclone® Anti-B (ABO2) and/or Seraclone® Anti-A,B (ABO3)

Materials required but not provided

- Pipettes (drop volume 40 to 50 µl)
- Isotonic saline solution
- Negative Control (e.g. Biotest Seraclone® Control ABO+Rh [REF] 805171100)
- Glass tubes 10 x 75mm or 12 x 75mm
- Serological Centrifuge
- Interval Timer
- Markers
- Optical aid (optional). The use of an optical aid for agglutination reading must be validated by the user.

Test Procedure

Tube test

1. Prepare a 3 to 5% suspension of red blood cells to be tested in isotonic saline.
2. Place one drop reagent into an appropriately labeled tube.
3. Add one drop of red blood cell suspension into the tube labeled for it and mix.
4. Centrifuge for 20 seconds at 800 -1000 x g.
5. Gently dislodge red blood cell button and observe for agglutination.
6. Record results

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

