

Blood Grouping Reagent

Anti- \bar{s} (MNS4)

Seraclone[®] Human Monoclonal
(P3YAN3)

FOR IN-VITRO DIAGNOSTIC USE

For Tube Testing

MEETS FDA POTENCY REQUIREMENTS

U.S. License Number: 1798

Package size

[REF] 808068100 [VOL] 2 mL Seraclone[®] Anti- \bar{s} (MNS4)

Intended Use

For the determination of the \bar{s} (MNS4) antigen of red blood cells using the tube test.

Summary

Antibodies to the \bar{s} antigen usually occur following immunization and are capable of causing hemolytic disease of the fetus and newborn (HDFN) and hemolytic transfusion reactions (HTR)¹. The complex system of the MNS system consists of over 40 antigens carried on two glycoprotein molecules. M, N, S, \bar{s} , and U antigens are the most important antigens of the MNS system with regard to transfusion medicine.

The frequencies of the common phenotypes are shown in the table.

Phenotypes and Frequencies in the MNS System ¹		
Phenotype	Whites	Blacks
M+N-	28	26
M+N+	50	44
M-M+	22	30
S+s-U+	11	3
S+s+U+	44	28
S-s+U+	45	69
S-s-U-	0	Less than 1
S-s-U+w	0	Rare*

* May not be detected by some reagent and are listed as U-

Biotest Anti-M, Anti-N, Anti-S, and Anti- \bar{s} Blood Group Reagents are used to test for the presence or absence of the M, N, S, and \bar{s} antigens. They are used principally in the resolution of antibody problems or in family studies.

Principle of the Test

The test principle is hemagglutination. The antibody in Seraclone[®] Anti- \bar{s} (MNS4) binds to the antigen on red blood cells. This does not result in a direct agglutination reaction. By adding Anti-Human Globulin reagent the antibody coated red blood cells are linked to each other, visible as red blood cell agglutination.

Reagent

As the reactive component Seraclone[®] Anti- \bar{s} (MNS4) contains a human monoclonal antibody of the immunoglobulin class IgG. It is derived from cell culture supernatant and demonstrates the consistent specificity and red blood reproducibility characteristic for monoclonal antibodies.

Antibodies are diluted in a buffered protein solution containing macromolecular potentiators.

The following antibody is produced using intermediate products produced for Biotest Medical Diagnostics GmbH in a shared manufacturing agreement with DIAGAST, Parc Eurasante, 251 av. Eugene Avinee-BP9, 59374 Loos Cedex France; License Number 1744.

Seraclone[®] Anti- \bar{s} (MNS4) clone P3YAN3 (IgG)

Preservative: 0.1% sodium azide, 0.05g/L sodium arsenite

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 all 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- \bar{s} (MNS4)

Materials required but not provided

- Pipettes (drop volume 40 to 50 μ l)
- Isotonic saline solution
- Anti-Human Globulin Anti-IgG (e.g. Biotest [REF] 804175100)
- IgG coated red blood cells (e.g. Biotest Coombscell-E [REF] 816030100)
- Glass tubes 10 x 75mm or 12 x 75mm
- Serological Centrifuge
- IgG coated red blood cells
- 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 and mix.
4. Incubate for 30 minutes at 37°C.
5. Wash red blood cells three times with isotonic saline solution. Completely decant the supernatant.
6. Follow the directions of the Anti-Human Globulin manufacturer.
7. Centrifuge for 20 seconds at 800 -1000 x g.
8. Gently dislodge red blood cell button and observe for agglutination.
9. Record results

Stability of the Reaction

Following centrifugation, all tube tests should be read immediately and results interpreted without delay. Time delays may cause a dissociation of the antigen-antibody complexes resulting to false negative or more often weak positive reactions.

Quality Control

The reactivity of all blood typing reagents should be confirmed by testing with known positive and negative red blood cells on each day of use.

To confirm the reactivity or specificity of Biotest Monoclonal Anti- \bar{s} Blood Grouping Reagent, it should be tested with antigen-positive (preferably from heterozygous individuals) and antigen-negative red blood cells, respectively. The reagent is satisfactory for use if it reacts only with antigen-positive red blood cells.

Negative results in an antiglobulin test should be verified with IgG coated red blood cells: Add one drop of IgG coated red blood cells, mix and centrifuge for 20 seconds at 800 - 1000 x g. Positive result: The

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

negative reaction in the indirect antiglobulin test is valid, reactive Anti-Human Globulin is present. Negative result: A technical error was made and the test must be repeated.

It is recommended that a positive and a negative control be performed in parallel with testing.

Interpretation of results

Agglutination of the red blood cells is a positive result and indicates the presence of the corresponding antigen. No agglutination is a negative result and indicates the absence of the corresponding antigen.

An agglutination viewer may facilitate the reading of tube tests (as recommended by the AABB Technical Manual, 15th edition).

Frequencies in the population are listed in the "Summary" section.

Limitations

- Samples with a positive direct antiglobulin test, cold agglutinins, or rouleaux formation may show false positive results in testing with monoclonal antibodies. Results on these samples must be interpreted with caution. False positive results or reaction suspected to be due to cold agglutinins should be resolved according to in-house procedures. It is recommended that an appropriate control be tested in parallel.
- Insufficient or inappropriate washing can lead to false negative or false positive reactions. Small amounts of residual patient sera/plasma can neutralize the Anti-Human Globulin.
- Some conditions that may cause false positive results are:
 - Contamination of sample or reagents
 - Autoantibodies
 - Improper storage or preparation of red blood cells
 - Antibodies to antibiotics or other reagents
 - Cold Antibodies

Specific Performance Characteristics

Testing is performed in accordance with FDA recommended methods. The final release testing is performed according to the product specific SOPs. Each lot of Biotest Blood Group Reagent is tested in the Quality control by package insert method against a panel of antigen positive red blood cells (heterozygous antigen expression and if possible weakened antigen expression) to insure suitable reactivity. The products meet FDA potency requirements. The specificity testing for the presence of contaminating antibodies is performed according to the product specific SOPs.

For the product performance it is necessary to adhere to the recommended method in the instructions for use.

The performance of the Biotest Anti- \bar{s} was confirmed against a FDA approved reference reagent in a Multi Center Field Trial.

For Technical Support or further product information, contact Biotest Diagnostics Corporation at 800-522-0090.

Note

Each facility should verify the optimum spin time for the specific centrifuge in use.

Manual techniques are to be performed according to the manufacturer's instructions. Each deviation from these instructions is the sole responsibility of the user.

Used tests must be discarded as hazardous material. Manage waste according to local, state and national regulations.

Glossary of Symbols

Symbol	Definition	Symbol	Definition
LOT	Batch Code	IVD	<i>In vitro</i> diagnostic medical device
⚠	Caution, consult accompanying documents	📖	Consult instructions for use.
🏭	Manufacturer	🕒	Use by YYYY-MM-DD
⚠	Contains sufficient quantity for <n> tests.	REF	Catalog number
🌡	Temperature limitation	VOL	Volume

Bibliography

1. Mark E. Brecher, MD et al. Technical Manual 15th Edition, Bethesda, MA: AABB, 2005.
2. Marion E. Reid, Christine Lomas-Francis, The Blood Group Antigen FactsBook, New York, NY: Academic Press, 2004.

For Reference Use Only