

**Culture supernatant Mouse Anti-human  
Blood Group A Antibody \*HE-195,  
monoclonal\***Catalog number: V103260  
Unit size: 1.0 ml**Product Details**

---

|                    |                                                                     |
|--------------------|---------------------------------------------------------------------|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze.                 |
| Expiration Date    | 12 months upon receiving                                            |
| Concentration      | Lot specific (please consult certificate of analysis for given lot) |
| Formulation        | Culture supernatant, 0.2% (w/v) BSA                                 |

**Antibody Properties**

---

|                    |                     |
|--------------------|---------------------|
| Species Reactivity | Human               |
| Class              | Primary             |
| Clonality          | Monoclonal          |
| Host               | Mouse               |
| Immunogen          | Blood Group A       |
| Clone              | HE-195              |
| Conjugate          | Culture supernatant |

**Biological Properties**

---

|             |                                                                                                                    |
|-------------|--------------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with Culture supernatant under optimal conditions |
| Application | AGG                                                                                                                |

**Applications**

---

Identifies erythrocytes and glycoproteins of blood group A.