

**PE Mouse Anti-human IgG (Fab) Antibody
*4A11, monoclonal***

Catalog number: V1031480

Unit size: 0.1 mg

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 | Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA |

Antibody Properties

| | |
|--------------------|------------|
| Species Reactivity | Human |
| Class | Secondary |
| Clonality | Monoclonal |
| Host | Mouse |
| Immunogen | IgG (Fab) |
| Clone | 4A11 |
| Conjugate | PE |

Biological Properties

| | |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions |
| Application | FC (QC TESTED) |

Applications

Detects the antigen-binding fragment of Immunoglobulin G.